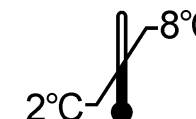
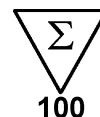













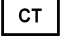

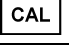

ALDOSTERONE



ALDO-RIACT



<p>Trousse pour le dosage radioimmunologique de l'aldostérone dans le sérum, le plasma ou l'urine</p> <p>Pour diagnostic In Vitro</p> <p>La trousse contient :</p> <table border="0"> <tr> <td>Tubes revêtus</td> <td>2 x 50 tubes</td> </tr> <tr> <td>Traceur ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Calibrateur 0</td> <td>1 x qsp 4 mL</td> </tr> <tr> <td>Calibrateurs 1 – 5</td> <td>5 x qsp 1 mL</td> </tr> <tr> <td>Contrôle</td> <td>1 x qsp 1 mL</td> </tr> <tr> <td>Notice d'utilisation</td> <td>1</td> </tr> </table> <p>Attention : Certains réactifs contiennent de l'azoture de sodium</p>	Tubes revêtus	2 x 50 tubes	Traceur ≤ 111 kBq	1 x 50 mL	Calibrateur 0	1 x qsp 4 mL	Calibrateurs 1 – 5	5 x qsp 1 mL	Contrôle	1 x qsp 1 mL	Notice d'utilisation	1	<p>Kit for the radioimmunoassay of aldosterone in serum, plasma or urine</p> <p>For In Vitro diagnostic use</p> <p>Kit content :</p> <table border="0"> <tr> <td>Coated tubes</td> <td>2 x 50 tubes</td> </tr> <tr> <td>Tracer ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Calibrator 0</td> <td>1 x qs 4 mL</td> </tr> <tr> <td>Calibrators 1 – 5</td> <td>5 x qs 1 mL</td> </tr> <tr> <td>Control</td> <td>1 x qs 1 mL</td> </tr> <tr> <td>Instruction for use</td> <td>1</td> </tr> </table> <p>Warning : Some reagents contain sodium azide</p>	Coated tubes	2 x 50 tubes	Tracer ≤ 111 kBq	1 x 50 mL	Calibrator 0	1 x qs 4 mL	Calibrators 1 – 5	5 x qs 1 mL	Control	1 x qs 1 mL	Instruction for use	1	<p>Kit zur radioimmunologischen Bestimmung von Aldosteron in Serum, Plasma oder Urin</p> <p>Zur In Vitro Diagnostik</p> <p>Inhalt des Kits :</p> <table border="0"> <tr> <td>Teströhrchen beschichtet</td> <td>2x 50 Röhrchen</td> </tr> <tr> <td>Tracer ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Kalibrator 0</td> <td>1 x qs 4 mL</td> </tr> <tr> <td>Kalibratoren 1 – 5</td> <td>5 x qs 1 mL</td> </tr> <tr> <td>Kontrolle</td> <td>1 x qs 1 mL</td> </tr> <tr> <td>Gebrauchsinformation</td> <td>1</td> </tr> </table> <p>Achtung : Einige Reagenzien enthalten Natriumazid</p>	Teströhrchen beschichtet	2x 50 Röhrchen	Tracer ≤ 111 kBq	1 x 50 mL	Kalibrator 0	1 x qs 4 mL	Kalibratoren 1 – 5	5 x qs 1 mL	Kontrolle	1 x qs 1 mL	Gebrauchsinformation	1
Tubes revêtus	2 x 50 tubes																																					
Traceur ≤ 111 kBq	1 x 50 mL																																					
Calibrateur 0	1 x qsp 4 mL																																					
Calibrateurs 1 – 5	5 x qsp 1 mL																																					
Contrôle	1 x qsp 1 mL																																					
Notice d'utilisation	1																																					
Coated tubes	2 x 50 tubes																																					
Tracer ≤ 111 kBq	1 x 50 mL																																					
Calibrator 0	1 x qs 4 mL																																					
Calibrators 1 – 5	5 x qs 1 mL																																					
Control	1 x qs 1 mL																																					
Instruction for use	1																																					
Teströhrchen beschichtet	2x 50 Röhrchen																																					
Tracer ≤ 111 kBq	1 x 50 mL																																					
Kalibrator 0	1 x qs 4 mL																																					
Kalibratoren 1 – 5	5 x qs 1 mL																																					
Kontrolle	1 x qs 1 mL																																					
Gebrauchsinformation	1																																					
<p>kit per il dosaggio radioimmunologico dell'aldosterone nel siero, nel plasma o nell'urina</p> <p>Per uso diagnostico In Vitro</p> <p>Contenuto del kit :</p> <table border="0"> <tr> <td>Provette coattate</td> <td>2 x 50 provette</td> </tr> <tr> <td>Tracciante ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Calibratore 0</td> <td>1 x q.b a 4 mL</td> </tr> <tr> <td>Calibratori 1 – 5</td> <td>5 x q.b a 1 mL</td> </tr> <tr> <td>Controllo</td> <td>1 x q.b a 1 mL</td> </tr> <tr> <td>Istruzioni per l'uso</td> <td>1</td> </tr> </table> <p>Attenzione : Alcuni reagenti contengono sodio azide</p>	Provette coattate	2 x 50 provette	Tracciante ≤ 111 kBq	1 x 50 mL	Calibratore 0	1 x q.b a 4 mL	Calibratori 1 – 5	5 x q.b a 1 mL	Controllo	1 x q.b a 1 mL	Istruzioni per l'uso	1	<p>Equipo para la determinación inmunnoradiométrica cuantitativa específica de aldosterona en suero, plasma u orina</p> <p>Para uso diagnóstico In Vitro</p> <p>Contenido del equipo :</p> <table border="0"> <tr> <td>Tubos recubiertos</td> <td>2 x 50 tubos</td> </tr> <tr> <td>Trazador ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Calibrador 0</td> <td>1 x csp 4 mL</td> </tr> <tr> <td>calibradores 1 – 5</td> <td>5 x csp 1 mL</td> </tr> <tr> <td>Control</td> <td>1 x csp 1 mL</td> </tr> <tr> <td>Instrucciones de uso</td> <td>1</td> </tr> </table> <p>Precauciones : Algunos reactivos contienen azida sódica</p>	Tubos recubiertos	2 x 50 tubos	Trazador ≤ 111 kBq	1 x 50 mL	Calibrador 0	1 x csp 4 mL	calibradores 1 – 5	5 x csp 1 mL	Control	1 x csp 1 mL	Instrucciones de uso	1	<p>Immunoradiometrisk analysats för aldosteron i serum, plasma eller urin</p> <p>För invitro diagnostik</p> <p>Innehåll:</p> <table border="0"> <tr> <td>Belagda rör</td> <td>2 x 50 rör</td> </tr> <tr> <td>Spårämnet ≤ 111 kBq</td> <td>1 x 50 mL</td> </tr> <tr> <td>Kalibratur 0</td> <td>1 x 4 mL</td> </tr> <tr> <td>Kalibraturer 1 – 5</td> <td>5 x 1 mL</td> </tr> <tr> <td>Kontroll</td> <td>1 x 1 mL</td> </tr> <tr> <td>Bruksanvisning</td> <td>1</td> </tr> </table> <p>Varning: vissa reagenser innehåller natriumazid</p>	Belagda rör	2 x 50 rör	Spårämnet ≤ 111 kBq	1 x 50 mL	Kalibratur 0	1 x 4 mL	Kalibraturer 1 – 5	5 x 1 mL	Kontroll	1 x 1 mL	Bruksanvisning	1
Provette coattate	2 x 50 provette																																					
Tracciante ≤ 111 kBq	1 x 50 mL																																					
Calibratore 0	1 x q.b a 4 mL																																					
Calibratori 1 – 5	5 x q.b a 1 mL																																					
Controllo	1 x q.b a 1 mL																																					
Istruzioni per l'uso	1																																					
Tubos recubiertos	2 x 50 tubos																																					
Trazador ≤ 111 kBq	1 x 50 mL																																					
Calibrador 0	1 x csp 4 mL																																					
calibradores 1 – 5	5 x csp 1 mL																																					
Control	1 x csp 1 mL																																					
Instrucciones de uso	1																																					
Belagda rör	2 x 50 rör																																					
Spårämnet ≤ 111 kBq	1 x 50 mL																																					
Kalibratur 0	1 x 4 mL																																					
Kalibraturer 1 – 5	5 x 1 mL																																					
Kontroll	1 x 1 mL																																					
Bruksanvisning	1																																					

	FRA	ENG	DEU	ITA	SPA	POR	SWE	POL	HUN	CES	BUL	RUS	SRB	ELL
	Explication des symboles	Explanation of symbols	Erläuterung der Symbole	Spiegazione dei simboli	Significado de los símbolos	Explicação dos símbolos	Symbol förklaring	Wyjaśnieni e symboli	Jelmagyarázat	Vysvětlení symbolů	Объяснение на символите	Объяснение на символите	Objašnjenje simbola	Επεξήγηση συμβόλων
	Conforme aux normes européennes	European conformity	CE-Konformitätskennzeichnung	Conformità europea	Conformidad europea	conformidade Europeia	Förenlig med europeiska normer	Zgodne z normami europejskimi	Megfelel az európai szabványoknak	Evropská shody	Европейската съответствие	Европейский соответствия	Evropska usaglašenost	Επεξήγηση συμβόλων
	T° limite de stockage	Storage temperature limitation	Limitierung der Lagertemperatur	Limiti per la temperatura di conservazione	Limites de temperatura de almacenamiento	Limitação de temperatura de armazenamente	T°-gräns vid förvaring	Graniczna temperatura przechowywania	Tárolási hőmérséklet határ	Mezní teplota skladování	Ограничаване на температурата на съхранение	Ограничение температуры хранения	Ograničenje temperature za čuvanje	Περιορισμός θερμοκρασίας αποθήκευσης
	N° de lot	Batch code	Chargencode	codice lotto	Código de lote	código do lote	Lotnr.	Numer partii	Gyártási szám	Č. šarže	номер	код партии	Šifra serije	Κωδικός партиδας
	Utiliser jusqu'au	Use by	Verwendbar bis	utilizzare entro	Consumir antes de	usado por	Används senast	Zużyć do	Felhasználható az alábbi dátumig :	Použitelné do	Срок на валидност	Используйте по	Upotrebiti do	Χρήση με
	Consulter la notice d'utilisation	Consult operating instructions	Das Handbuch zu Rate ziehen	consultare le istruzioni per l'USO	Consultar las instrucciones de manejo o funcionamiento	Consulte o manual de instruções	Läs bruksanvisningen	Patrz dołączona ulotka	Olvassa el a használati utasítást	Přečtěte si návod k použití	Консултирайте инструкциите за работа	Обратитесь к инструкции по эксплуатации	Pogledajte uputstvo za upotrebu	Συμβουλευτείτε τις οδηγίες λειτουργίας
	Diagnostic In Vitro	In Vitro Diagnostic device	In-Vitro Diagnostische Anwendung	Dispositivo Diagnostico In Vitro	Dispositivo de diagnóstico In Vitro	No dispositivo de diagnóstico in vitro	In vitro-diagnos	Diagnostyka In Vitro	In vitro diagnosztika	Diagnostika in vitro	За ин витро диагностика устройство	Экстракорпоральное диагностическое устройство	Uređaj za dijagnostiku <i>in vitro</i>	Διαγνωστική συσκευή In Vitro
	Fabriqué par	Manufactured by	Hergestellt von	Prodotto da	Fabricado por	fabricado por	Tillverkad av	Wyprodukowane przez	Gyártja:	Vyrobil	Произведено от	Изготовитель	Proizveo	Κατασκευάζεται από
	Référence	Catalogue number	Katalog Nr.	N. catalogo	Número de catálogo	número de catálogo	Referens	Wzorzec	Referenciakészítmény	Reference	Каталожен номер	номер по каталогу	Kataloški broj	Αριθμός καταλόγου
	Nombre de tubes	Number of determinations	Anzahl der Bestimmungen	Numero di determinazioni	Número de determinaciones	Número de determinações	Antal rör	Liczba próbek	A kémcsövek száma	Počet zkumavek	Брой определяния	Количество определений	Broj određivanja	Αριθμός προσδιορισμών
	Tubes revêtus	Coated tubes	beschichtete Röhrchen	Provette coattate	Tubos recubiertos	tubos revestidos	Belagda rör	Probówki powlekane	Bevont kémcsövek	Zkumavky	Покритите тръби	Покрытые трубы	Obložene epruvete	Επικαλυμμένοι σωλήνες
	Traceur radioactif	Radioactive tracer	Radioactiver Tracer	Traccianti radioattivo	Trazador radiactivo	traçador radioativo	Radioaktiv tracer	Znacznik radioaktywny	Nyomjelző izotóp	Tracer	Индикатор	радиоактивного индикатора	Radioaktivni indikator	Ραδιενεργός ιχνηλάτης
	Calibrateur	Calibrator	Kalibrator	Calibratore	Calibrador	Calibrador	Kalibrator	Kalibrator	Kalibrátor	Kalibrátor	Калибратор	калибратор	Kalibrator	Πρότυπο
	Contrôle	Control	Kontrolle	Controllo	Control	controle	Kontroll	Kontrola	Kontroll	Kontrola	Контрол	контроль	Kontrola	έλεγχος

FRA

Modifications par rapport à la version précédente :
Information interférence biotine, ajout du Grec.

ENG

Changes from the previous version:
Biotin interference information, addition of Greek.

DEU

Änderungen gegenüber der Vorgängerversion:
Biotin Interferenzinformation, Zugabe von Griechisch.

ITA

Modifiche rispetto alla versione precedente:
Informazioni sull'interferenza della biotina, aggiunta di Greco.

SPA

Cambios desde la versión anterior:
Información de interferencia de biotina, Además de Griego.

POR

Alterações em relação à versão anterior:
Informações sobre interferência de biotina, adição de Grego.

SWE

Ändringar från föregående utgåva:
Biotininterferensinformation, tillägg av Grekiska.

POL

Zmiany w stosunku do poprzedniej wersji:
Informacja o zakłóceniach biotyiny, dodatek Grecki.

HUN

Változások az előző verzióhoz képest:
Biotin interferencia információ, Görög hozzáadásával.

CES

Změny od předchozí verze:
Informace o interferenci s biotinem, doplněk řečtiny.

BUL

Промени от предишната версия :
Информация за смущения в биотина, добавяне на гръцки.

RUS

Изменения по сравнению с предыдущей версией:
Информация об интерференции биотина, добавление греческого

SRB

Izmene u odnosu na prethodnu verziju:
Informacije o interferenciji biotina, dodavanje Grčkog.

ELL

Αλλαγές από την προηγούμενη έκδοση:
Πληροφορίες παρεμβολής βιοτίνης, προσθήκη ελληνικών.

1. NAME AND INTENDED USE

ALDO-RIACT is a kit for the radioimmunoassay of aldosterone in serum, plasma or urine. The kit is intended for professional use.

2. INTRODUCTION

Aldosterone is a steroid hormone with a molecular weight of 360.4. It is secreted by the glomerulosa zone of the adrenal gland and controls regulation of hydromineral metabolism.

In the distal part of the nephron, aldosterone encourages reabsorption of Na⁺ and Cl⁻ ions and secretion of K⁺ and H⁺ ions in blood vessel lumen. Thus there is an increase in extra-cellular osmolarity and water retention.

Regulation of aldosterone secretion is linked to the body's water balance and brings three factors into play. These are the renin-angiotensin system, the plasmic Na⁺/K⁺ ratio and, as a secondary factor, the ACTH.

Physiologically, there seems to be a circadian rhythm for levels of aldosterone in the blood, which may be linked to changes in body posture. It has also been found that aldosterone levels are higher for children than for adults, and increase during pregnancy.

In pathology, lowered levels are found in cases of adrenal insufficiency and in some congenital adrenal hyperplasia cases with enzyme deficiencies. However, aldosterone assay is mainly indicated for complete etiological examinations related to arterial hypertension (AHT), when a high level of aldosterone is a sign of hyperaldosteronism, either primary (adrenal adenoma) or secondary (malignant AHT, renal artery stenosis...).

3. PRINCIPLE

The principle of the assay is based on competition between a fixed amount of 125 iodine aldosterone and the aldosterone contained in calibrators or specimens to be assayed for a fixed and limited number of antibody binding sites bound to the solid phase (coated tubes).

After incubation and removal of the unbound tracer, the amount of 125 iodine aldosterone bound to the antibody is inversely proportional to the amount of unlabelled aldosterone present in the sample.

4. REAGENTS

Each kit contains enough reagents for 100 tubes. The expiry date is marked on the external label.

REAGENTS	SYMBOLS	QUANTITY	STORAGE
COATED TUBES : ready for use Anti-aldosterone antibody coated on the bottom of the tube.	CT	2 packs of 50 tubes	2-8°C until the expiry date.
ALDOSTERONE ¹²⁵I : ready for use ¹²⁵ I labelled aldosterone, buffer, bovine albumin, rabbit serum, sodium azide, red dye, non specific immunoglobulins. ≤ 111 kBq (≤ 3 μCi) per vial	TRACER	1 50 mL vial	2-8°C until the expiry date.
CALIBRATOR 0 : lyophilized Human aldosterone-free serum, preservative. Reconstitute the vial's contents with 4 mL of distilled water.	CAL	1 qs 4 mL vial	2-8°C until the expiry date. After reconstitution: 1 week at 2-8°C or 1 month at -20°C within the limits of the expiry date of the kit.
CALIBRATORS : lyophilized** Human serum, preservative, aldosterone 25, 60, 180, 500, 1500 pg/mL* (69, 166, 498, 1385, 4150 pmol/L). Reconstitute the vial's contents with 1 mL of distilled water.	CAL	5 qs 1 mL vials	2-8°C until the expiry date. After reconstitution: 1 month at - 20°C within the limits of the expiry date of the kit.
CONTROL : lyophilized** Human serum, preservative, aldosterone 110 pg/mL* (304 pmol/L). Reconstitute the vial's contents with 1 mL of distilled water.	CONTROL	1 qs 1 mL vial	2-8°C until the expiry date. After reconstitution: 1 month at - 20°C within the limits of the expiry date of the kit.

(*) The values shown above are only target values: the true value of each calibrator or control is shown on its label.

(**) Calibrators and control should be frozen and thawed only once.

5. PRECAUTIONS FOR USE

5.1. Safety measures

Raw materials of human origin contained in the reagents of this kit have been tested with licensed kits and found negative for the anti-HIV 1, anti-HIV 2, anti-HCV antibodies and the HBs antigen. However as it is impossible to strictly guarantee that such products will not transmit hepatitis, the HIV virus, or any other viral infection, all raw materials of human origin including the samples to be assayed must be treated as potentially infectious.

Do not pipette by mouth.

Do not smoke, eat or drink in areas in which specimens or kit reagents are handled.

Wear disposable gloves while handling kit reagents or specimens and wash hands thoroughly afterwards.

Avoid splashing.

Decontaminate and dispose of specimens and all potentially contaminated materials as if they contained infectious agents. The recommended method of doing this is autoclaving for a minimum of one hour at 121.5°C.

Sodium azide may react with lead or copper piping to form highly explosive metal azides. During waste disposal, flush the drains thoroughly to prevent a build-up of these products.

5.2. Basic radioprotection rules

This radioactive product may only be received, purchased, stored or used by persons so authorized, and by laboratories covered by such authorization. The solution should under no circumstances be administered to humans or to animals.

The purchase, storage, use or exchange of radioactive products are subject to the laws in force in the user's country.

Enforcement of the basic radioprotection rules will ensure adequate security.

A summary of these is given below:

Radioactive products must be stored in their original containers in a suitable area.

A record of the reception and storage of radioactive products must be kept up to date.

Handling of radioactive products should take place in a suitably-equipped area with restricted access (controlled zone).

Do not eat, drink, smoke or apply cosmetics in a controlled zone.

Do not mouth-pipette radioactive solutions.

Avoid any direct contact with all radioactive products by using laboratory coats and protective gloves.

Contaminated laboratory equipment and glassware must be disposed of immediately after contamination to prevent cross-contamination of different isotopes.

Any contamination or radioactive substance loss should be dealt with in accordance with the established procedures.

All radioactive waste disposal must be carried out according to the regulations in force.

5.3. Handling precautions

Do not use kit components beyond their expiry date.

Do not mix reagents from different batches.

Do not manipulate much than 100 tubes at the same time.

Avoid any microbic contamination of the reagents or of the water.

Fully respect the incubation times and the washing instructions indicated.

6. SPECIMEN COLLECTION AND PREPARATION

6.1. Serum or plasma

The assay is performed directly on serum or plasma; haemolyzed or hyperlipemic samples should not be used. If the test is to be carried out within 24 hours, the samples must be refrigerated at 2-8°C. Otherwise, they should be divided into aliquots and deep frozen (-20°C) until needed.

Dilutions

Should elevated aldosterone levels be suspected, the Calibrator 0 found in the kit is used for dilution.

It is recommended to carry out the dilutions using disposable plastic tubes.

6.2. Urine

Collect 24-hour urine, measure and record the volume. Mix well before drawing off an aliquot to assay. Add 1g boric acid / 100 mL of urine and store at 2-8 °C, or at - 20°C for extended storage.

Urine hydrolysis

Aldosterone can be assayed in urine samples after acid hydrolysis of aldosterone 18-glucuronide. In the conditions recommended for the assay, hydrolysis is complete.

- Mix 500 µL urine and 1 mL 0.1 N HCl. Use glass tubes.

- Cap tubes and incubate for 15-20 hours at 30 ± 2°C.

7. ASSAY PROCEDURE

7.1. Material required

Precision micropipettes or similar, with disposable tips, capable of dispensing 10 µL, 200 µL, 500 µL, 1 mL and 4 mL. Their calibration should be checked regularly.

Distilled water.

Disposable plastic tubes.

Vortex-type mixer.

Circular horizontal shaker (400 rpm).

Gamma scintillation counter calibrated for 125 iodine measurement.

7.2. Protocol

All reagents must be brought to room temperature (18-25°C) at least 30 minutes before their use. Calibrators and control must be reconstituted 15 minutes before use. Dispensing of the reagents into the tubes is carried out at room temperature (18-25°C).

The assay requires the following groups of tubes:

Calibrator "0" group, for the determination of maximum binding (Bo).

Calibrator groups, to establish the calibrator curve.

Control group for the control.

Sx groups, to test serum or plasma samples.

Ux groups, to test urine samples.

It is recommended that the assay be performed in duplicate for the calibrators, the control and the samples.

Strictly respect the order in which reagents are to be added:

Calibrators, control or serum and plasma samples

Dispense 200 µL of calibrators, control or samples to be assayed into the corresponding tubes.

Urine samples

Dispense 200 μ L of calibrator 0 + 10 μ L of hydrolysed urine samples to be assayed into the corresponding tubes.

Add 0.5 mL of 125 I Aldosterone to each tube.

Mix each tube gently with a Vortex-type mixer.

Incubate for 3 h \pm 5 mn at room temperature (18-25°C) under shaking (400 rpm).

Aspirate the contents of the tubes as completely as possible. Any trace of red dye should disappear.

Measure the remaining radioactivity bound to the tubes with a gamma scintillation counter calibrated for 125 iodine measurement.

8. QUALITY CONTROL

Good laboratory practices require that quality control samples be used in each series of assays to check the quality of the results obtained. All specimens should be treated identically, and result analysis using the appropriate statistical methods is recommended.

9. RESULTS

For each group of tubes, calculate the mean counts after subtracting the background.

Calculate the B/Bo ratio for each calibrator and unknown sample as follows:

$$B/Bo (\%) = \frac{\text{CAL cpm or Sx cpm or Ux cpm}}{\text{CAL 0 cpm}} \times 100$$

Draw up the calibrator curve by plotting the calibrators' B/Bo against their concentrations.

Serum or plasma

Read the sample values directly from the curve, correcting the read value for the dilution factor, if necessary.

Urine

Multiply the read value on the calibration curve by 63 to obtain the urinary aldosterone concentration in pg/mL. The factor 63 results from the dilution factor of hydrolyzed urine (factor 3) multiplied by the (calibrator+sample)-to-sample ratio (factor 21).

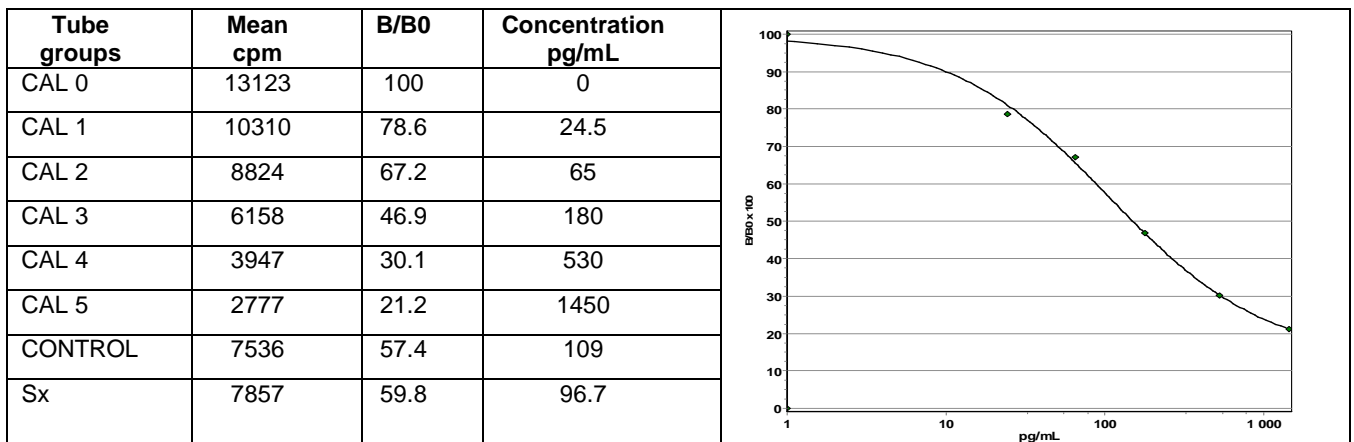
Daily aldosterone excretion expressed as μ g of aldosterone/24 hours is calculated thus:

$$\text{pg/mL} \times \text{mL urine/24 h} \times 10^{-6} = \mu\text{g/24 hours}$$

Conversion of μ g aldosterone in nmol is: 1 μ g = 2.77 nmol.

The "forced" spline mathematical fitting model is recommended for calibration curve. Other fitting model may give slightly different results.

Typical calibrator curve (example only): this data must not be substituted for results obtained in the laboratory.



10. PROCEDURAL LIMITATIONS

Samples containing fibrin, gross haemolysis, gross lipemia or turbidity may give misleading results.

Falsely elevated levels may be obtained for patients undergoing substitute corticotherapy.

Do not attempt to extrapolate sample values beyond the last calibrator. Dilute the samples and retest.

11. EXPECTED VALUES

11.1. Plasma expected values:

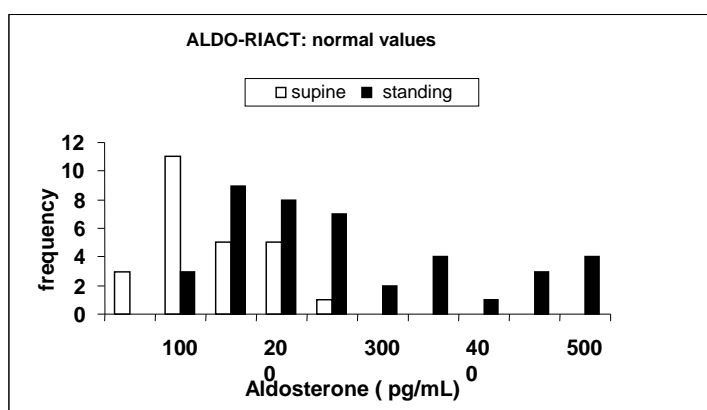
The ranges given below are only indicative for a normal population with a normal sodium intake; each laboratory should establish its own reference ranges.

NORMAL SUBJECTS	PLASMA (pg/mL)		
	5 th percentile	Median	95 th percentile
Supine position* (n=25)	42	99	201.5
Standing position (n=41)	97	201	626

* In supine position for at least one hour.

Conversion of pg/mL aldosterone in pmol/L is: 1 pg/mL = 2.77 pmol/L.

Because of possible serie interferences, this method without preliminary sample extraction can give some different values compared to an extraction method.



11.2. Urinary expected values:

The ranges given below are only indicative for a normal population with a normal sodium intake; each laboratory should establish its own reference ranges.

NORMAL SUBJECTS	URINARY VALUES (µg / 24 h)			
	Mean	Standard deviation	Minimum	Maximum
n = 110	4.7	3.75	0.12	17.6

12. SPECIFIC CHARACTERISTICS OF THE ASSAY

12.1. Imprecision

This has been assessed using 2 samples with different concentrations. They were tested either 29 times in the same series of assays, or in duplicate in 15 different series.

Sample	Mean pg/ml	Within-run CV %	Between-run CV %
1	30.1	7.7	8.4
2	656	8.3	5.0

12.2. Specificity

The antibody used in this assay guarantees a measurement which is completely specific for aldosterone. It shows very low cross-reactions with some other compounds.

Aldosterone	100 %	Dexamethasone	N.D. %
Androstene-dione	0.00217 %	Estradiol	0.00037 %
Androsterone	0.00065 %	Estriol	N.D. %
Dehydroepi-Androsterone	0.00016 %	Estrone	0.00019 %
Canrenone	N.D. %	Prazosin-HCl	N.D. %
Cortexolone (11-deoxycortisol)	0.0052 %	Prednisolone	0.00035 %
Corticosterone	0.04 %	Prednisone	0.00022 %
11-deoxy-Corticosterone	0.065 %	Pregnanetriol	N.D. %
18-hydroxydeoxy-Corticosterone	0.012 %	Pregnenolone	0.00043 %
Cortisol	0.0022 %	Progesterone	0.027 %
Cortisone	0.0018 %	17 α -hydroxy-Progesterone	0.0025 %
9 α -fludro-Cortisone	N.D. %	Spironolactone	0.00037 %
		Testosterone	0.0026 %

N.D. = Not detectable

12.3. Detection limit

The detection limit is defined as being the smallest detectable concentration different from zero with a probability of 95 %. It has been assessed as being 7 pg/mL.

12.4. Measuring range

7 – 1500 pg/mL

12.5. Interferences

- Biotin interference was measured from 1.5 ng/mL.
- No interference with haemoglobin and triglycerides, measured up to respective concentrations of equal to 2.5 g/L and 10 g/L has been observed.

The **ALDO-RIACT** kit is protected against heterophilic antibodies. However, we cannot guarantee that this cover is complete.

ASSAY FLOW CHART

Tubes	Calibrator 0 μ L	Calibrators Control Serum or Plasma Samples μ L	Hydrolysed urine samples μ L	Aldosterone ¹²⁵ I μ L	Mix Incubate 3 H \pm 5 mn at 18-25°C under shaking Aspirate	Count
Calibrator 0	200			500		
Calibrators		200		500		
Control		200		500		
Serum Sx		200		500		
Urine Ux	200		10	500		

Bibliography

Balikian HM, Brodie AH, Dale SL, Melby JC, Tait JF. Effect of posture on the metabolic clearance rate, plasma concentration and blood production rate of aldosterone in man. *J Clin Endocrinol Metab.* 1968;28:1630-40.

Conn JW. Presidential address: Part I. Painting background. Part II. Primary aldosteronism, a new clinical syndrome. *J Lab Clin Med.* 1955;45:3-17.

Ganguly A, Dowdy AJ, Luetscher JA, Melada GA. Anomalous postural response of plasma aldosterone concentration in patients with aldosterone-producing adrenal adenoma. *J Clin Endocrinol Metab.* 1973;36:401-4.

Grimaux M, Fernandez B, Malard V, Seguin P. Monoclonal antibody-based radioimmunoassay of aldosterone. Proceedings of the Ninth International Congress of Endocrinology. Nice: International society of endocrinology; 1992:P04.31.032.

Hegstad R, Brown RD, Jiang NS, et al. Aging and aldosterone. *Am J Med.* 1983;74:442-8.

Nomura K, Han DC, Jibiki K, Demura H, Tsushima T, Shizume K. Primary aldosteronism with normal aldosterone levels in blood and urine. *Acta Endocrinol.* 1985;110:522-5.

Schambelan M, Brust NL, Chang BCF, Slater KL, Biglieri EG. Circadian rhythm and effect of posture on plasma aldosterone concentration in primary aldosteronism. *J Clin Endocrinol Metab.* 1976;43:115-31.

Young WF, Hogan MJ. Primary aldosteronism: Diagnosis and treatment. *Mayo Clin Proc.* 1990;65:96-110