



Thermo Scientific Clinical Chemistry Reagent Values For The Pointe Scientific Chemistry Calibrator





REF C7506-50

LOT 832502


 2011-12

Cal. **LOT** 832502  2012-10

Dil. **LOT** 832502  2011-12



Constituent / Constituant Bestandteil / Costituente Componente / Συστατικό	Method / Méthode / Methode Metodo / Método / Μέθοδος			Value / Valeur / Wert Valore / Valor / Αξία	
				(SI) 37°C	(non SI) 37°C
ALB* Albumin / Albumine / Albumina / Albúmina / ευκωματίνη	BCG	TR360		36 g/L	3.6 g/dL
BICARB* Bicarbonate / Bikarbonat Bicarbonato / διπτανθρακικό	Enzymatic / Enzymatique / Enzymatische/ Enzimatico Enzimático / Ενζυματικός	Infinity™ TR283		18 mmol/L	18 mEq/L
		Infinity™ TR282		17 mmol/L	17 mEq/L
BILI-T* Total Bilirubin Bilirubine Totale Gesamtbilirubin Bilirubina Totale Bilirubina Total Ολικής Χολερυθρίνης	Surfactant / Agent de surface Tensid / Surfattanti Tensioactivo Επιφανειοδραστικά	TR323		70 µmol/L	4.1 mg/dL
		1245		83 µmol/L	4.8 mg/dL
				72 µmol/L	4.2 mg/dL
BILI-D** Direct Bilirubin / Bilirubine Directe Direktbilirubin / Bilirubina Diretta Bilirubina Directa άμεσης Χολερυθρίνης	Diazo	TR333		49 µmol/L	2.9 mg/dL
CA* Calcium Calcio Ασβεστίου	Arsenazo III Αρσενάζο III	TR292		2.55 mmol/L	10.2 mg/dL
		Infinity™ 1265		2.58 mmol/L	10.3 mg/dL
CHOL* Cholesterol / Cholestérol Cholesterin / Colesterolo Colesterol / Χοληστερόλη	Enzymatic CHO/CHE / Enzymatique CHO/CHE Enzymatische CHO/CHE / Enzimatico CHO/CHE Enzimático CHO/CHE / Ενζυματικός CHO/CHE	TR133		5.04 mmol/L	195 mg/dL
		Infinity™ TR135			
		Infinity™ TR134			
CRE* Creatinine Créatinine Kreatinin Creatinina κρεατινίνη	Kinetic Jaffe / Cinétique Jaffe Kinetisch Jaffe / Cinetico Jaffe Cinético Jaffe Κινητική Jaffe	Infinity™ TR351		472 µmol/L	5.34 mg/dL
		Enzymatic / Enzymatique / Enzymatische Enzimatico / Enzimático / Ενζυματικός	TR354		375 µmol/L
GLU* Glucose Glukose Glucosio Glucosa Γλυκόζη	Hexokinase Esochinasi Hexoquinasa Εξοκινάση	TR150		10.11 mmol/L	182 mg/dL
		Infinity™ TR154		10.39 mmol/L	187 mg/dL
		Infinity™ TR152		10.07 mmol/L	181 mg/dL
		TR151		10.51 mmol/L	189 mg/dL

© 2010 Thermo Fisher Scientific Inc. All rights reserved.

 Fisher Diagnostics
a division of Fisher Scientific Company, LLC
a part of Thermo Fisher Scientific Inc.
Middletown, VA 22645-1905 USA
Phone: 800-528-0494
540-869-3200
Fax: 540-869-8132

REF C7506-50

10 x 5 mL
1 x 55 mL

Constituent / Constituant Bestandteil / Costituente Componente / Συστατικό	Method / Méthode / Methode Metodo / Método / Μέθοδος			Value / Valeur / Wert Valore / Valor / Αξία	
				(SI) 37°C	(non SI) 37°C
FE* Iron / Fer / Eisen Ferro / Hierro / Σιδήρου	Ferrozine	TR461		29 µmol/L	165 µg/dL
Mg* Magnesium Magnésium Magnesio Μαγνήσιο	Arsenazo / Αρσενάζο	Infinity™ TR311		1.51 mmol/L	3.03 mEq/L
PHOS** Inorganic Phosphorus Phosphore Inorganique Anorganischer Phosphor Fosforo Inorganico Fósforo Inorgánico Ανόργανος Φώσφορος	UV/Ammonium molybdate UV/Molybdate d'ammonium UV/Molibdato di ammonio UV/Molibdato de amonia UV/Ammonium-molybdat UV/Μολυβδαινικό αμμώνιο	TR300		1.81 mmol/L	5.6 mg/dL
TP* Total Protein / Protéines Totales Gesamtprotein / Proteine Totali Πρωτεΐνες Totales / Ολική Πρωτεΐνη	Biuret	TR340		60 g/L	6.0 g/dL
TG* Triglycerides Triglycérides Triglyzeride Trigliceridi Triglicéridos Τριγλυκερίδια	GPO	TR222		1.16 mmol/L	102 mg/dL
		Infinity™ TR223			
		Infinity™ TR224		1.04 mmol/L	92 mg/dL
UREA NITROGEN* Urea Urée Harnstoff Ουρία	UV-Kinetic / UV-Cinétique UV-Kinetisch / UV-Cinetico UV-Cinético UV-Κινητική	TR120		16.5 mmol/L	46 mg/dL
		Infinity™ TR123		16.5 mmol/L	46 mg/dL
		Infinity™ TR124			
UA* Uric Acid Acide Urique Harnsäure Acido Urico Ácido Úrico Ουρικό Οξύ	Enzymatic/Trinder Enzymatique/Trinder Enzymatische/Trinder Enzimatico/Trinder Enzimático/Trinder Ενζυματικός/Trinder	TR240		426 µmol/L	7.2 mg/dL
		Infinity™ TR242		430 µmol/L	7.2 mg/dL
		Infinity™ TR243			

*Known concentrations verified against available NIST Standard Reference Material.

** Value is not NIST verified.



For in vitro diagnostic use. Utilisation en diagnostique in vitro. Für in vitro Diagnostik. Per uso diagnostico in vitro. Para uso en diagnósticos in vitro. Προοριζόμενο για διάγνωση in vitro.



Batch code/Lot number. Numéro de lot. Batch Code/Losnummer. Codice/Numero lotto. Código de lote/Número de lote. Αριθμός παρτίδας.



Catalogue number. Référence catalogue. Katalognummer. Numero catalogo. Número de catálogo. Αριθμός Καταλόγου.



Consult instructions for use. Consulter les instructions d'utilisation. Siehe Benutzungsvorschriften. Consultare le istruzioni d'uso. Consulte las instrucciones de uso. Συμβουλευτείτε τις οδηγίες χρήσης.



Temperature Limitation. Limites de température. Temperaturbeschränkung. Limite di temperatura. Limitación de temperatura. Περιορισμός θερμοκρασίας.



Use by/Expiration Date. Utiliser jusque. Verfallsdatum. Usare entro/Data di scadenza. Usar hasta/Fecha de caducidad. Χρήση μέχρι/ημερ/νια λήξης.



Calibrator Reference Serum. Calibreur Sérum de Référence. Kalibrator Referenzserum. Calibratore Siero di Riferimento. Calibrador Suero de Referencia. Πρότυπος Ορός Βαθμονομητής.



CAUTION. CONSULT INSTRUCTIONS FOR USE. ATTENTION: Consulter les instructions d'utilisation. VORSICHT. Siehe Benutzungsvorschriften. AVVERTENZA. Consultare le istruzioni d'uso. PRECAUCIÓN. Consulte las instrucciones de uso. ΠΡΟΣΟΧΗ: Συμβουλευτείτε τις οδηγίες χρήσης.



Manufactured by. Fabriqué par. Hergestellt von. Prodotto da. Fabricado por. Κατασκευασμένο από.



With sample blanking. Avec test blanc de l'échantillon. Mit Probenleerwert. Con blanking del campione. Con blanco de muestra. Με τύφλωση του δείγματος.



Without sample blanking. Sans test blanc de l'échantillon. Without Probenleerwert. Senza blanking del campione. Sin blanco de muestra. Χωρίς τύφλωση του δείγματος.



Monochromatic. Monochromatique. Monochromatisch. Monocromatico. Monocromático. Μονοχρωματικός.



Bichromatic. Bichromatique. Bichromatisch. Bicromatico. Bicromático. Διχρωματικός.