









# Data-Cal™ Chemistry Calibrator

REF TS106500

LOT 430802-313  2017-12-31

Cal.  430802-313  2018-09-30

Dil.  430802-313  2017-12-31

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
<b>ALB*</b> Albumin / Albumine / Albumina / Albúmina / ευκωματίνης	BCG	TR360		37 g/L	3.7 g/dL
<b>BICARB*</b> Bicarbonate / Bikarbonat Bicarbonato / διπτανθρακικό	Enzymatic / Enzymatique / Enzymatische/ Enzimatico Enzimático / Ενζυματικός	Infinity™ TR283		18 mmol/L	18 mEq/L
		Infinity™ TR282		NA mmol/L	NA mEq/L
<b>BILI-T*</b> Total Bilirubin Bilirubine Totale Gesamtbilirubin Bilirubina Totale Bilirubina Total Ολικής Χολερυθρίνης	Van den Bergh (Liquid) / Van den Bergh (Liquide) Van den Bergh (Flüssig) / Van den Bergh (Liquido) Van den Bergh (líquido) / Van den Bergh (Υγρό)	1245		83 µmol/L	4.88 mg/dL
				71 µmol/L	4.18 mg/dL
<b>BILI-D**</b> Direct Bilirubin / Bilirubine Directe Direktbilirubin / Bilirubina Diretta Bilirubina Directa άμεσης Χολερυθρίνης	Diazo	TR333		45 µmol/L	2.65 mg/dL
<b>CA*</b> Calcium Calcio Ασβεστίου	Arsenazo III Αρσενάζιο III	Infinity™ 1265		2.42 mmol/L	9.7 mg/dL
<b>CHOL*</b> Cholesterol / Cholestérol Cholesterin / Colesterolo Colesterol / Χοληστερόλη	Enzymatic CHO/CHE / Enzymatique CHO/CHE Enzymatische CHO/CHE / Enzimatico CHO/CHE Enzimático CHO/CHE / Ενζυματικός CHO/CHE	Infinity™ TR134		4.63 mmol/L	179 mg/dL
<b>CRE*</b> Creatinine Creatinine Kreatinin Creatinina κρεατινίνης	Kinetic Jaffe / Cinétique Jaffe Kinetisch Jaffe / Cinetico Jaffe Cinético Jaffe Κινητική Jaffe	Infinity™ TR351		436 µmol/L	4.93 mg/dL
<b>GLU*</b> Glucose Glukose Glucosio Glucosa Γλυκόζη	Hexokinase Esocinasi Hexoquinasa Εξοκινάση	Infinity™ TR154		9.9 mmol/L	178 mg/dL
	Oxidase Οξειδάση	Infinity™ TR152		9.8 mmol/L	177 mg/dL
<b>FE*</b> Iron / Fer / Eisen Ferro / Hierro / Σιδήρου	Ferrozine	TR461		33 µmol/L	184 µg/dL
<b>TP*</b> Total Protein / Protéines Totales Gesamtprotein / Proteine Totali Proteínas Totales / Ολική Πρωτεΐνης	Biuret	TR340		59 g/L	5.9 g/dL
<b>TG*</b> Triglycerides Triglycérides Triglyzeride Trigliceridi Triglicéridos Τριγλυκερίδια	GPO	Infinity™ TR224		1.28 mmol/L	113 mg/dL

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
<b>UREA NITROGEN*</b> Urea Urée Harnstoff Ουρία	UV-Kinetic / UV-Cinétique UV-Kinetisch / UV-Cinetico UV-Cinético UV-Κινητική	Infinity™ TR124		15.6 mmol/L	44 mg/dL
<b>UA*</b> Uric Acid Acide Urique Harnsäure Acido Úrico Ácido Úrico Ουρικό Οξύ	Enzymatic/Trinder Enzymatique/Trinder Enzymatische/Trinder Enzimatico/Trinder Enzimático/Trinder Ενζυματικός/Trinder	Infinity™ TR243		367 µmol/L	6.2 mg/dL

\*Known concentrations verified against available NIST Standard Reference Material.

\*\* Value is not NIST verified.

© 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



MDCI Ltd.  
Arundel House  
1 Liverpool Gardens  
Worthing, West Sussex BN11 1SL UK



**Thermo**  
SCIENTIFIC