

In-vitro Diagnostic Reagents

N Terminal Procollagen 3 Peptide

Product number: N Terminal Procollagen 3 **Lot number:** 16B

Reagents	Peptide	Batch #	Expiry Date
Réactifs		Lot N°	Péremption
CONJ-P3NP-ELISA		16A	21/01/2025
MICROPLATE-P3NP-ELIS		16A	14/01/2025
DIL CAL0-P3NP-ELISA		16A	24/01/2025
CAL1-P3NP-ELISA		16A	29/01/2025
CAL2-P3NP-ELISA		16A	29/01/2025
CAL3-P3NP-ELISA		16A	29/01/2025
CAL4-P3NP-ELISA		16A	29/01/2025
CAL5-P3NP-ELISA		16A	29/01/2025
CONT1-P3NP-ELISA		16A	29/01/2025
CONT2-P3NP-ELISA		16A	29/01/2025
TWEEN 20		216A	18/06/2026
BLISTER-3-WASH		26A	31/03/2025
SUBS TMB		59AP	27/01/2025
STOP SOLN		35AP	07/02/2025
N Terminal Procollag		010	

Results / Résultats

Parameter	Conc.	Unit	Response	Measured value
Paramètre	Conc.	Unité	Réponse	Valeur mesurée
CAL 0	0.00	µg/L	DO	0.06
CAL 1	3.50	µg/L	DO	0.41
CAL 2	6.70	µg/L	DO	0.73
CAL 3	12.60	µg/L	DO	1.18
CAL 4	21.60	µg/L	DO	1.84
CAL 5	32.00	µg/L	DO	2.25

Controls	Unit	Expected values	Measured value
Contrôles	Unité	Valeurs attendues	Valeur mesurée
CONTROL	µg/L	5.920 - 8.880	7.38
CONTROL 2	µg/L	14.800 - 22.200	18.13

Control Date 17/07/2023

Date du contrôle

Quality control manager visa

Visa du responsable du contrôle qualité

Expiry date 14/01/2025

Date de péremption



DELEUZE MAZZOCCA Isabelle