

SHBG-RIACT

Protocole/Protocol : 30mn + 30mn

Lot N°/Batch # : 222A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-SHBG-RIACT	222A	12/01/2022
CT-SHBG-RIACT	19A	16/01/2023
BUF-SHBG-RIACT	20A	19/08/2022
WASH-SHBG-RIACT	15A	18/08/2022
CAL0-SHBG-RIACT	22A	21/02/2023
CAL1-SHBG-RIACT	23A	09/05/2023
CAL2-SHBG-RIACT	23A	09/05/2023
CAL3-SHBG-RIACT	23A	09/05/2023
CAL4-SHBG-RIACT	23A	09/05/2023
CAL5-SHBG-RIACT	23A	09/05/2023
CONT1-SHBG-RIACT	22A	21/02/2023
CONT2-SHBG-RIACT	22A	21/02/2023
SHBG-RIACT MO	015	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	nmol/L	B/T %	0.06
CAL.A	4.40	nmol/L	B/T %	3.41
CAL.B	16.60	nmol/L	B/T %	11.01
CAL.C	35.00	nmol/L	B/T %	23.14
CAL.D	80.00	nmol/L	B/T %	41.03
CAL.E	150.00	nmol/L	B/T %	56.00

Activité totale / Total activity : 90545 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	nmol/L	18.40 - 27.60	21.90
CONTROL 2	nmol/L	48.80 - 73.20	54.75

Date du contrôle
Control Date

15/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

12/01/2022

