

Batch information

Mab GST-XL

HTRF MAb Anti GST-XL665 - 20K pts

Product reference :

61GSTXLB

Product characteristics :

- Lot Number	39A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.12.09
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Mab)	400.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	400.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	20 000

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

