

Mab GST-XL

MAB Anti GST-XL665 - 20,000 tests

Product reference :

61GSTXLB**Product characteristics :**

- Lot Number_____	36A
- 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_____	2022.10.01
- Reconstitute each vial with_____	1 mL distilled H2O
- Active Moiety Concentration (Mab)_____	400 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

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