

anti GST-d2

MAb Anti GST-d2 - 4 mg

Product reference :

61GSTDAJ

Product characteristics :

- Lot Number_____	08A
- Packaging Unit_____	200 000 tests/vial
- Conditioning buffer_____	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance_____	Frozen form
- Storage conditions_____	<-16 °C
- Expiration date_____	2025.07.26
- Volume per vial_____	8 mL
- Active Moiety Concentration (GST)_____	500 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (GST)	4.00 mg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (GST)	20.00 ng
Number of tests/vial	200 000

*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 HTRF web site: www.htrf.com/products