

STD

Mouse TNF alpha kit - Standard

Product reference :

62MTNFACDA

Product characteristics :

- Lot Number_____ 05A
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 1% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.11.27

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	5 400.00 pg	N/A

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