

STD

Human IgG kit standard

Product reference :

62HFCCDA

Product characteristics :

- Lot Number_____ 16A
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 0.2% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.01.07

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	4 000.00 ng	N/A

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

