

STD

Human MMP9 kit - Standard
62MM9CDA

Product reference :

Product characteristics :

- Lot Number_____ 03A
- Conditioning buffer_____ Phosphate 50 mM, BSA 0,
2% + stabilizers
- Appearance_____ Frozen form
- Storage conditions_____ < -16 °C
- Expiration date_____ 2023.09.25
- Volume per vial_____ 50 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ag)	200.00 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