

**plasma sample diluent - 15ml**

Plasma sample diluent - 15 mL

Product reference :

**62DLPDDD**

**Product characteristics :**

- Lot Number	11A
- Conditioning buffer	-
- Appearance	Liquid form
- Storage conditions	4 °C
- Expiration date	2023.06.18
- Volume per vial	100 mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	15.00 mL	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](http://www.cisbio.com)*