

Batch information

STD

HTRF cAMP Gs HiRange kit - Std

Product reference :

62AM6CDA

Product characteristics :

- Lot Number	15A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 0.2% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.10.16

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (AMP)	6 720.00 pmoles	351

PLEASE NOTE

The expiration date of batch 15A has been extended to 2024-10-16 after validation by QC

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

