

Batch information

Buffer enz HTRF Enzymatic Buf./kinase (5X) - 10 mL
Product reference : **62EZBFDC**

Product characteristics :

- Lot Number 14A
- Conditioning buffer -
- Appearance Liquid form
- Storage conditions 2-8 °C
- Expiration date 2026.11.18

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	10.00 mL	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

