

Batch information

Control lysate HTRF BTK p-Y223 kit - Ctrl lysate
Product reference : **63ADK017TDA**

Product characteristics :

- Lot Number 06A
- Conditioning buffer Lysis Buffer
- Appearance Frozen form
- Storage conditions -60 °C
- Expiration date 2025.01.17
- Volume per vial 150.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	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

