

Batch information

control lysate

HTRF ATF2 p-T71 kit - Ctrl lysate

Product reference :

63ADK015TDA**Product characteristics :**

- Lot Number	06A
- Conditioning buffer	Phosphate 0.1M pH7 + % BSA + % Tween
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.10.12

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	150 000

PLEASE NOTE

The expiration date of batch 06A has been extended to 2024-10-12 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

