

## Batch information

**SAH d2 conjugate - Medium**

HTRF Methyltransferase kit - 10K pts d2

Product reference :

**62SAHDLA**

### Product characteristics :

- Lot Number	13A
- Conditioning buffer	H2O, 0.1% BSA
- Appearance	Liquid form
- Storage conditions	< -60 °C
- Expiration date	2025.03.24
- Volume per vial	5.50 mL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SAH)	5.50 mL	N/A

### PLEASE NOTE

See package insert of #62SAHPEH for use.

***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)***

