

D2V2-NHS

d2v2-NHS - 1 mg

Product reference :

65DVSABB

Product characteristics :

- Lot Number _____ 06A
- Conditioning buffer _____ ///

- Appearance _____ Desiccated form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2022.09.26

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (DV) | 1.00 mg | 1 066 |

PLEASE NOTE

Rz (D0650nm/280nm)=34.7

Reconstitute with DMSO

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

