

I-BET-762

Catalog #: 27630-1

Lot #: 141016

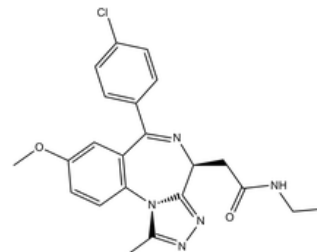
Size: 5 mg

Structure:

CAS Registry #: 1260907-17-2

Purity: ≥98%

Chemical Formula: C₂₂H₂₂ClN₅O₂



Molecular Weight: 423.9

Description: I-BET-762 is a small molecule inhibitor of the BET family of bromodomain-containing proteins. It has been shown to bind to the acetylated lysine recognition motifs on the bromodomain of BET proteins. This binding prevents the interaction of between BET proteins and acetylated histone peptides.

Appearance: White powder

Solubility: Soluble in DMSO at 85 mg/ml, water at 11 mg/ml, or ethanol at 85 mg/ml.

Biological Activity: I-BET-762 has a reported IC₅₀ of roughly 35 nM

Storage/Stability: Powder is stable at -20 °C for up to 3 years. Re-suspended powder is stable at -80 °C for up to 6 months.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Nicodeme, E., *et al.*, *Nature*. 2010. **468(7327)**: 1119-1123.
2. Bandukwala, H.S., *et al.*, *Proc. Nat. Acad. Sci.* 2012. **109(36)**: 14532-14537.