

Product Information

Description

Iberdomide is a derivative of Thalidomide and has antitumor and immunostimulatory activities because it is a potent cereblon (CRBN) E3 ligase modulator.

Molecular Formula:

$C_{25}H_{27}N_3O_5$

Supplied As:

Solid

Stability:

At least 6 months at -20°C . Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

449.5 Da

Cas No.

1323403-33-3

Chemical Structure

