

Product Information

Description

Pomalidomide is the third-generation immunomodulatory agent, functions through interacting with the E3 ligase cereblon and induces degradation of essential Ikaros transcription factors.

Molecular Formula:

C₁₃H₁₁N₃O₄

Supplied As:

Solid

Solubility:

DMSO

Stability:

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

Storage:

-20°C

MW:

273.2 Da

Cas No.

19171-19-8

Purity:

98%

Chemical Structure

