

Product Information

Description

Pomalidomide-PEG5-Azide is a PROTAC linker that incorporates the Pomalidomide based cereblon ligand and 4-unit PEG linker used in PROTAC technology. Azide enable click chemistry with Alkyne, DBCO or BCN molecules.

Molecular Formula:

$C_{26}H_{34}N_6O_{10}$

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

590.6 Da

Purity:

98%

Chemical Structure

