

Product Information

Description

3-(5-amino-2-methyl-4-oxoquinazolin-3(4H)-yl)piperidine-2,6-dione has been shown to treat Diffuse large B-cell lymphoma with antitumor and immunomodulatory activity. Its molecular target is the protein cereblon, a substrate receptor of the cullin ring E3 ubiquitin ligase complex CRL4.

Molecular Formula:

C₁₄H₁₄N₄O₃

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

286.3 Da

Cas No.

1015474-32-4

Chemical Structure

