

Product Information

Description

Tucatinib acts as an HER2 inhibitor, and in BT474 breast carcinoma cells, it effectively suppresses HER2 phosphorylation. Tucatinib inhibits AKT phosphorylation and induces apoptosis in BT474 cells, rendering it a viable cytotoxic agent for formulating antibody-drug conjugates (ADCs).

Molecular Formula:

$C_{26}H_{24}N_8O_2$

Supplied As:

Solid

Stability:

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

Storage:

-20°C

MW:

480.5 Da

Cas No.

937263-43-9

Chemical Structure

