

## Product Information

## Description

BETd-260 is a PROTAC connected by ligands for Cereblon and BET. It potently suppresses cell viability and induces apoptosis in hepatocellular carcinoma (HCC) cells.

## Molecular Formula:

$C_{43}H_{46}N_{10}O_6$

## Supplied As:

Solid

## Stability:

At least 6 months at  $-20^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

## Storage:

$-20^{\circ}\text{C}$

## MW:

798.9 Da

## Cas No.

2093388-62-4

## Chemical Structure

