

## Product Information

## Description

Thalidomide is a drug used in the treatment of erythema nodosum leprosum (ENL) and multiple myeloma. It can inhibit cereblon (CRBN) and be used in the context of targeted protein degradation as molecular glue or PROTAC.

## Molecular Formula:

$C_{13}H_{10}N_2O_4$

## Supplied As:

Solid

## Stability:

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

## Storage:

-20°C

## MW:

258.2 Da

## Cas No.

50-35-1

## Purity:

98%

## Chemical Structure

