

## Product Information

## Description

ARV-825 is a PROTAC molecule composed of a BRD4 binding moiety joined to an E3 ligase cereblon binding moiety. ARV-825 recruits BRD4 to E3 ubiquitin ligase cereblon, resulting in the degradation of BRD4.

## Molecular Formula:

$C_{46}H_{47}ClN_8O_9S$

## Supplied As:

Solid

## Stability:

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

## Storage:

$-20^{\circ}C$

## MW:

923.5 Da

## Cas No.

1818885-28-7

## Chemical Structure

