

## Product Information

### Description

I-[(3R)-3-[4-amino-3-(4-phenoxyphenyl)-1Hpyrazolo[3,4-d]pyrimidin-1-yl]-1-piperidinyl]-2-propyn-1-one, is a derivative of the first-in-class covalent BTK inhibitor ibrutinib. Ibrutinib inhibits B-cell proliferation and survival by irreversibly binding the protein Bruton's tyrosine kinase (BTK). B cell cancers are directly affected by this since the B-cell receptor pathway is often aberrantly active.

### Molecular Formula:

C<sub>25</sub>H<sub>22</sub>N<sub>6</sub>O<sub>2</sub>

### Supplied As:

Solid

### Stability:

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

### Storage:

-20°C

### MW:

438.5 Da

### Cas No.

1970122-88-3

## Chemical Structure

