

Product Information

Description	IBT6A is a Btk kinase inhibitor. This compound is a precursor for the manufacture of ibrutinib.
Molecular Formula:	C ₂₂ H ₂₂ N ₆ O
Supplied As:	Solid
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	386.5 Da
Cas No.	1022150-12-4
Purity:	97%

Chemical Structure

