

UNC-2170

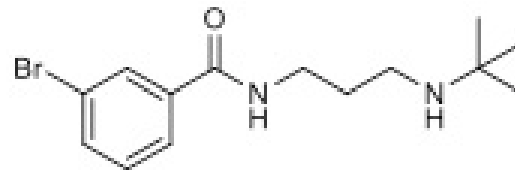
Catalog No: 14120, 14121**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 349.7

 $C_{14}H_{21}BrN_2O \cdot HCl$

CAS 1648707-58-7

Physical Properties: White powder; MP = 162-164°C

Names: UNC-2170, 3-Bromo-N-(3-(tert-butylamino)propyl)benzamide, hydrochloride

Pharmacology: The DNA damage response protein 53BP1 utilizes its tandem tudor domain to recognize dimethylated lysine 20 on histone (H4K20me2), a modification associated with double-stranded DNA breaks. UNC-2170, identified by screening, was found to be a μ M ligand for 53BP1 which also demonstrated at least 17-fold selectivity for 53BP1 over other methyl-lysine binding proteins. The compound functions as 53BP1 antagonist in cell lysates and suppresses class switch recombination in whole cells, a process requiring a functioning 53BP1 tudor domain (ref 1).

Solubilization: May be dissolved in DMSO (40 mg/ml); or water (50 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. MT Perfetti *et al.* *ACS Chem. Biol.* 2015, 10:1072

Storage and Guarantee: Store desiccated as supplied at -20° for up to 2 years. Store solutions at -20°C for up to 3 months. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.