

PRODUCT DATA SHEET

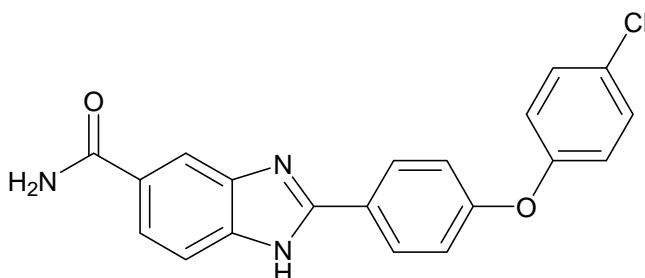


PRODUCT: BML-277 CAS NO.: 516480-79-8
2-(4-(4-chlorophenoxy)phenyl)-1H-benzimidazole-5-carboxamide
Chk2 inhibitor II

CATALOG NO.: EI-388

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Off-white solid

MOLECULAR FORMULA: $C_{20}H_{14}ClN_3O_2$ MOLECULAR WEIGHT: 363.8

PURITY: 98% (TLC: 50% Ethyl acetate/hexanes; $R_f = 0.17$)

SOLUBILITY: Soluble in: DMSO (20 mg/mL); ethanol (20 mg/mL)

STORAGE: Store, as supplied, at $-20^{\circ}C$ for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: BML-277 is a highly selective inhibitor of Chk2 ($IC_{50}=15$ nM) displaying <25% inhibition of 35 other kinases at $10 \mu M$ ¹. It displays potent radioprotective activity and prevents apoptosis of human T-cells subjected to ionizing radiation ($EC_{50}=3-7.6 \mu M$)¹. It is a useful tool for dissecting the role of Chk2 in cellular signaling².

REFERENCES: 1. K.L. Arienti *et al.* *J. Med. Chem.* 2005 **48** 1873
2. Y. Pereg *et al.* *Mol. Cell. Biol.* 2006 **26** 6819

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.