

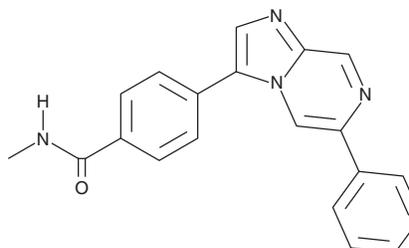
# PRODUCT INFORMATION



## AZ 32

Item No. 26380

**CAS Registry No.:** 2288709-96-4  
**Formal Name:** N-methyl-4-(6-phenylimidazo[1,2-a]pyrazin-3-yl)-benzamide  
**MF:** C<sub>20</sub>H<sub>16</sub>N<sub>4</sub>O  
**FW:** 328.4  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

AZ 32 is supplied as a solid. A stock solution may be made by dissolving the AZ 32 in the solvent of choice, which should be purged with an inert gas. AZ 32 is soluble in the organic solvent DMSO at a concentration of approximately 150 mg/ml.

### Description

AZ 32 is an inhibitor of ataxia-telangiectasia mutated (ATM) kinase (IC<sub>50</sub> = <6.2 nM).<sup>1</sup> It is selective for ATM over PI3Kα and ataxia-telangiectasia and Rad3-related (ATR) kinase (IC<sub>50</sub>s = 4.6 and >4.6 μM, respectively). AZ 32 enhances radiation-induced cytotoxicity in a panel of five human glioma cells expressing wild-type or mutant p53. It increases survival in a U87-281G glioma orthotopic mouse xenograft model, as well as a NCI H2228 non-small cell lung cancer (NSCLC) mouse xenograft model of metastatic brain tumors, when administered at doses of 200 and 50 mg/kg, respectively, in combination with ionizing radiation.

### Reference

1. Karlin, J., Allen, J., Ahmad, S.F., *et al.* Orally bioavailable and blood-brain barrier-penetrating ATM inhibitor (AZ32) radiosensitizes intracranial gliomas in mice. *Mol. Cancer Ther.* **17**(8), 1637-1647 (2018).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/23/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM