

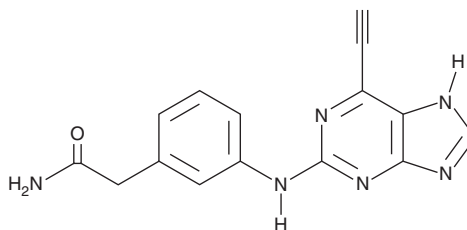
PRODUCT INFORMATION



NCL-00017509

Item No. 31462

CAS Registry No.: 1507367-00-1
Formal Name: 3-[[6-ethynyl-9H-purin-2-yl]amino]-benzeneacetamide
MF: $C_{15}H_{12}N_6O$
FW: 292.3
Purity: $\geq 98\%$
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

NCL-00017509 is supplied as a solid. A stock solution may be made by dissolving the NCL-00017509 in the solvent of choice, which should be purged with an inert gas. NCL-00017509 is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

Description

NCL-00017509 is an inhibitor of the serine/threonine kinase Nek2 ($IC_{50} = 56 \text{ nM}$).^{1,2}

References

1. Carbain, B., Bayliss, R., Boxall, K., *et al.* 2-Arylamino-6-ethynylpurines as potent irreversible inhibitors of the mitotic kinase Nek2. *Poster Session - Combinatorial Chemistry, Drug Delivery, Drug Design and Drug Synthesis* **118** (2012).
2. Lebraud, H., Coxon, C.R., Archard, V.S., *et al.* Model system for irreversible inhibition of Nek2: Thiol addition to ethynylpurines and related substituted heterocycles. *Org. Biomol. Chem.* **12**(1), 141-148 (2014).

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, 12/08/2022

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