

PRODUCT INFORMATION



GCN2-IN-1

Item No. 36027

CAS Registry No.: 1448693-69-3

Formal Name: 3-(1H-indazol-6-yl)-N-[1-(tetrahydro-2H-pyran-4-yl)-1H-pyrazol-4-yl]-3H-1,2,3-triazolo[4,5-d]pyrimidin-5-amine

Synonyms: A-92, GCN2 Inhibitor 1, General Control Nonderepressible 2 Kinase Inhibitor 1

MF: $C_{19}H_{18}N_{10}O$

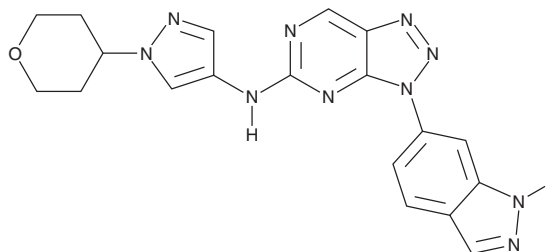
FW: 402.4

Purity: $\geq 90\%$

Supplied as: A crystalline 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

GCN2-IN-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the GCN2-IN-1 in the solvent of choice, which should be purged with an inert gas. GCN2-IN-1 is soluble in DMSO.

Description

GCN2-IN-1 is an inhibitor of general control nonderepressible 2 kinase (GCN2; $IC_{50} = \sim 0.2 \mu\text{M}$ for autophosphorylation), a serine/threonine protein kinase involved in the integrated stress response.¹ It also inhibits integrase phosphorylation ($IC_{50} = \sim 0.2 \mu\text{M}$ for the HIV-1 enzyme). GCN2-IN-1 inhibits ELAV-like protein 1, also known as human antigen R (HuR), protein dimerization ($IC_{50} = 4.5 \mu\text{M}$).² GCN2-IN-1 also decreases the viability of U251, XD456, U87, and LN-229 glioma cells (IC_{50} s = 4.7, 2.8, 3.2, and $2.7 \mu\text{M}$, respectively).

References

1. Torres, C., Garling, A., Taouji, S., *et al.* Targeting the integrated stress response kinase GCN2 to modulate retroviral integration. *Molecules* **26**(17), 5423 (2021).
2. Filippova, N., Yang, X., Ananthan, S., *et al.* Targeting the HuR oncogenic role with a new class of cytoplasmic dimerization inhibitors. *Cancer Res.* **81**(8), 2220-2233 (2021).

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, 10/04/2022

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