

# PRODUCT INFORMATION



**GS-444217**

Item No. 40757

**CAS Registry No.:** 1262041-49-5

**Formal Name:** 4-(4-cyclopropyl-1H-imidazol-1-yl)-N-[3-(4-cyclopropyl-4H-1,2,4-triazol-3-yl)phenyl]-2-pyridinecarboxamide

**Synonyms:** Apoptosis Signal-regulating Kinase 1-IN-1, Apoptosis Signal-regulating Kinase 1 Inhibitor 1, ASK1-IN-1, ASK1 Inhibitor 1

**MF:** C<sub>23</sub>H<sub>21</sub>N<sub>7</sub>O

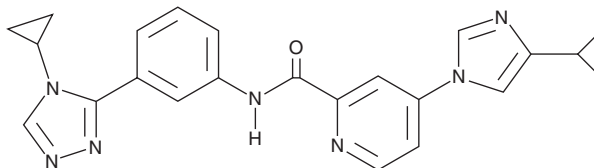
**FW:** 411.5

**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

GS-444217 is supplied as a solid. A stock solution may be made by dissolving the GS-444217 in the solvent of choice, which should be purged with an inert gas. GS-444217 is slightly soluble (0.1-1 mg/ml) in DMSO.

## Description

GS-444217 is an inhibitor of apoptosis signal-regulating kinase 1 (ASK1; IC<sub>50</sub> = 2.87 nM).<sup>1</sup> It is selective for ASK1 over 442 additional kinases at 1 μM. GS-444217 (0.001-10 μM) inhibits phosphorylation of the ASK1 substrates MAP kinase kinase 3 (MKK3), MKK6, and MKK4, as well as their downstream substrates p38 MAPK and JNK in HEK293T cells expressing human ASK1. Dietary administration of GS-444217 (0.3% w/w) reduces renal fibrosis and proteinuria in a *db/db eNOS<sup>-/-</sup>* mouse model of diabetic kidney disease (DKD).

## Reference

1. Liles, J.T., Corkey, B.K., Notte, G.T., *et al.* ASK1 contributes to fibrosis and dysfunction in models of kidney disease. *J. Clin. Invest.* **128**(10), 4485-4500 (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, 04/24/2024

## 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](http://WWW.CAYMANCHEM.COM)