

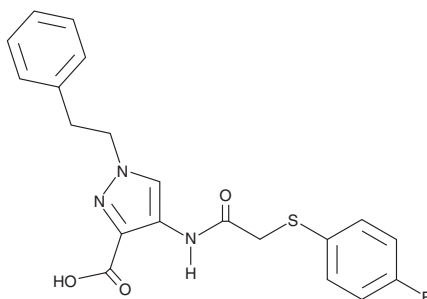
# PRODUCT INFORMATION



**KR-33493**

Item No. 22071

**CAS Registry No.:** 1021497-97-1  
**Formal Name:** 4-[[2-[(4-bromophenyl)thio]acetyl]amino]-1-(2-phenylethyl)-1H-pyrazole-3-carboxylic acid  
**MF:** C<sub>20</sub>H<sub>18</sub>BrN<sub>3</sub>O<sub>3</sub>S  
**FW:** 460.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 256 nm  
**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

KR-33493 is supplied as a crystalline solid. A stock solution may be made by dissolving the KR-33493 in the solvent of choice, which should be purged with an inert gas. KR-33493 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

## Description

KR-33493 is an inhibitor of Fas-associated factor 1 (FAF1).<sup>1,2</sup> It inhibits chemical hypoxia-induced apoptosis in H9c2 cardiomyocytes and hydrogen peroxide-induced cell death in SH-SY5Y cells when used at a concentrations of 10 and 20 μM, respectively.<sup>2,3</sup> KR-33493 (20 mg/kg per day) reduces decreases in striatal dopamine transporter activity in a mouse model of Parkinson's disease induced by MPTP.<sup>1</sup>

## References

1. Park, H.S., Moon, B.S., Yoo, S.-E., *et al.* Assessment of potential protective effect on dopaminergic neuron by an Fas-associated factor 1 inhibitor treatment using [<sup>18</sup>F]FE-PE2I PET in a Parkinson's disease mouse model. *J. Nucl. Med.* **58(Suppl 1)**, 13 (2017).
2. Yoo, S.-e., Yu, C., Jung, S., *et al.* Design and synthesis of fluorescent and biotin tagged probes for the study of molecular actions of FAF1 inhibitor. *Bioorg. Med. Chem. Lett.* **26(4)**, 1169-1172 (2016).
3. Jung, Y.-S., Kim, E., Kim, N., *et al.* Aminopyrazole derivatives, process for the preparation thereof, and composition for preventing or treating ischemic diseases containing the same. *Korea Research Institute of Chemical Technology. WO 2008/051047 A1*, (2008).

### 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, 11/07/2022

## 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