

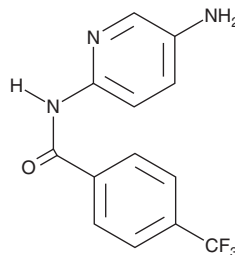
PRODUCT INFORMATION



TFAP

Item No. 26745

CAS Registry No.: 1011244-68-0
Formal Name: N-(5-amino-2-pyridinyl)-4-(trifluoromethyl)-benzamide
MF: $C_{13}H_{10}F_3N_3O$
FW: 281.2
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 223, 317 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

TFAP is supplied as a crystalline solid. A stock solution may be made by dissolving the TFAP in the solvent of choice, which should be purged with an inert gas. TFAP is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TFAP in DMF is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and ethanol.

TFAP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TFAP should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. TFAP has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TFAP is an inhibitor of COX-1 ($IC_{50} = 0.8 \mu\text{M}$).¹ It is selective for COX-1 over COX-2 ($IC_{50} = 210 \mu\text{M}$). TFAP (10 and 30 mg/kg) reduces acetic acid-induced writhing in mice and has analgesic activity in the formalin test in rats when administered at a dose of 30 mg/kg. It does not induce gastric damage in rats when administered at doses up to 300 mg/kg.

Reference

1. Kakuta, H., Zheng, X., Oda, H., *et al.* Cyclooxygenase-1-selective inhibitors are attractive candidates for analgesics that do not cause gastric damage. Design and in vitro/in vivo evaluation of a benzamide-type cyclooxygenase-1 selective inhibitor. *J. Med. Chem.* **51**(9), 2400-2411 (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, 10/24/2022

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