

PRODUCT INFORMATION



NSC 44625

Item No. 40985

CAS Registry No.: 654-62-6

Formal Name: 4-amino-6-(trifluoromethyl)-1,3-benzenedisulfonamide

Synonyms: Bendroflumethiazide Impurity A, 2,4-Disulfamyl-5-trifluoromethylaniline

MF: C₇H₈F₃N₃O₄S₂

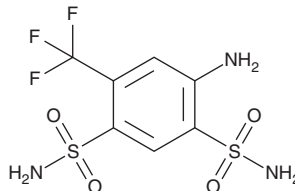
FW: 319.3

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

NSC 44625 is supplied as a solid. A stock solution may be made by dissolving the NSC 44625 in the solvent of choice, which should be purged with an inert gas. NSC 44625 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

NSC 44625 is an inhibitor of carbonic anhydrase II (CAII; K_i = 63 nM for the human enzyme).¹ It is selective for CAII over CAI (K_i = 5,800 nM for the human enzyme) but also inhibits β-carbonic anhydrase (K_i = 223 nM for the *S. cerevisiae* enzyme).

Reference

1. Isik, S., Kockar, F., Aydin, M., *et al.* Carbonic anhydrase inhibitors: inhibition of the β-class enzyme from the yeast *Saccharomyces cerevisiae* with sulfonamides and sulfamates. *Bioorg. Med. Chem.* **17**(3), 1158-1163 (2009).

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, 07/03/2024

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