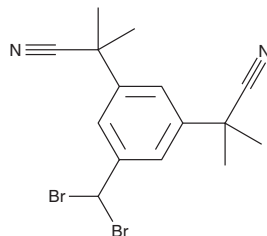


PRODUCT INFORMATION

5-Dibromomethyl Anastrozole

Item No. 36024

CAS Registry No.: 1027160-12-8
 Formal Name: 5-(dibromomethyl)- $\alpha^1, \alpha^1, \alpha^3, \alpha^3$ -tetramethyl-1,3-benzenediacetonitrile
 MF: $C_{15}H_{16}Br_2N_2$
 FW: 384.1
 Purity: $\geq 95\%$
 Supplied as: A solid
 Storage: $-20^\circ C$
 Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

5-Dibromomethyl anastrozole is supplied as a solid. A stock solution may be made by dissolving the 5-dibromomethyl anastrozole in the solvent of choice, which should be purged with an inert gas. 5-Dibromomethyl anastrozole is slightly soluble in chloroform and methanol.

Description

5-Dibromomethyl anastrozole is a synthetic intermediate in the synthesis of the aromatase/cytochrome P450 (CYP) isoform CYP19A1 inhibitor anastrozole (Item No. 11987).¹

Reference

1. Solanki, K.S., Pal, G., Haider, H., *et al.* Process for the preparation of pure anastrozole. **US2009/0221837A1** (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, 04/26/2023

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