

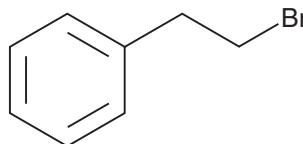
# PRODUCT INFORMATION



## Phenethylbromide

Item No. 20087

**CAS Registry No.:** 103-63-9  
**Formal Name:** (2-bromoethyl)-benzene  
**Synonym:** NSC 33926  
**MF:** C<sub>8</sub>H<sub>9</sub>Br  
**FW:** 185.1  
**Purity:** ≥98%  
**Supplied as:** A 10 mg/ml solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Phenethylbromide is an analytical reference material that is structurally categorized as an organobromide. It is used in the synthesis of a variety of compounds, including fentanyl. Phenethylbromide is combined with 4-piperidinone to produce N-phenethyl-4-piperidone, which, as a precursor in the synthesis of fentanyl, is scheduled as a List I chemical in the United States.<sup>1</sup> Phenethylbromide may be found as an impurity in samples of fentanyl produced using this pathway.

### Reference

1. D. E. A. D. o. Justice Control of a chemical precursor used in the illicit manufacture of fentanyl as a List I chemical. Final rule. *Federal Register* **73(144)**, 43355-43357 (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, 11/05/2025

#### 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