

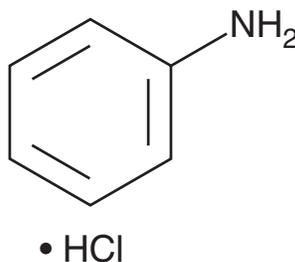
# PRODUCT INFORMATION



## Aniline (hydrochloride)

Item No. 21963

**CAS Registry No.:** 142-04-1  
**Formal Name:** benzenamine, monohydrochloride  
**Synonyms:** CI-76001, Phenylamine  
**MF:** C<sub>6</sub>H<sub>7</sub>N • HCl  
**FW:** 129.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 235, 288 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Aniline (hydrochloride) (Item No. 21963) is an analytical reference standard that is structurally categorized as a synthetic intermediate. It is used to make the imine derivative of NPP (Item No. 20528) as part of the synthesis of fentanyl (Item Nos. ISO60197 | 14719) and its derivatives.<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. Varadi, A., Palmer, T.C., Haselton, N., *et al.* Synthesis of carfentanil amide opioids using the ugi multicomponent reaction. *ACS Chem. Neurosci.* **6(9)**, 1570-1577 (2015).

#### 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, 06/13/2023

#### 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