

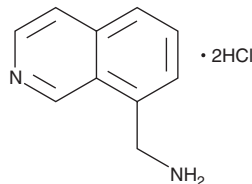
# PRODUCT INFORMATION



## 8-Isoquinoline methanamine (hydrochloride)

Item No. 10007125

**CAS Registry No.:** 850734-84-8  
**Formal Name:** 8-Isoquinolinemethanamine, dihydrochloride  
**MF:** C<sub>10</sub>H<sub>10</sub>N<sub>2</sub> • 2HCl  
**FW:** 231.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217 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

8-Isoquinoline methanamine (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of 8-isoquinoline methanamine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 8-isoquinoline methanamine (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

8-Isoquinoline methanamine (hydrochloride) is a synthetic intermediate useful for pharmaceutical synthesis.

#### 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, 10/13/2022

#### 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