

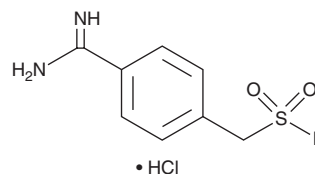
# PRODUCT INFORMATION



## *p*-APMSF (hydrochloride)

Item No. 14971

**CAS Registry No.:** 74938-88-8  
**Formal Name:** 4-(aminoiminomethyl)-benzenesulfonyl fluoride, monohydrochloride  
**Synonyms:** (*p*-Amidinophenyl)methanesulfonyl Fluoride, *para*-APMSF  
**MF:** C<sub>8</sub>H<sub>9</sub>FN<sub>2</sub>O<sub>2</sub>S • HCl  
**FW:** 252.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 232 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

*p*-APMSF (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the *p*-APMSF (hydrochloride) in the solvent of choice, which should be purged with an inert gas. *p*-APMSF (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of *p*-APMSF (hydrochloride) in DMSO and water is approximately 50 and 25 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

### Description

*p*-APMSF is an irreversible inhibitor of serine proteases with K<sub>i</sub> values of 1.02, 1.18, 1.5, and 1.54 μM for bovine trypsin, human thrombin, bovine plasmin, and bovine Factor Xa, respectively.<sup>1</sup> It is selective for these proteases over bovine chymotrypsin and acetylcholinesterase.

### Reference

1. Laura, R., Robison, D.J., and Bing, D.H. (*p*-Amidinophenyl)methanesulfonyl fluoride, an irreversible inhibitor of serine proteases. *Biochemistry* **19**(21), 4859-4864 (1980).

#### 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, 11/01/2022

#### 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