

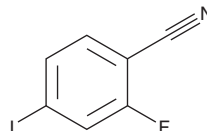
PRODUCT INFORMATION



2-Fluoro-4-iodo benzonitrile

Item No. 30483

CAS Registry No.: 137553-42-5
Synonym: 4-Iodo-2-fluoro benzonitrile
MF: C₇H₃FIN
FW: 247.0
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 280, 289 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-Fluoro-4-iodo benzonitrile is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-fluoro-4-iodo benzonitrile in the solvent of choice, which should be purged with an inert gas. 2-Fluoro-4-iodo benzonitrile is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-fluoro-4-iodo benzonitrile in these solvents is approximately 30 mg/ml.

2-Fluoro-4-iodo benzonitrile is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-fluoro-4-iodo benzonitrile should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 2-Fluoro-4-iodo benzonitrile has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2-Fluoro-4-iodo benzonitrile is a building block.^{1,2} It has been used in the synthesis of *L. infantum* trypanothione reductase (Li-TryR) dimerization and oxidoreductase activity inhibitors.¹ 2-Fluoro-4-iodo benzonitrile has also been used in the synthesis of transient receptor potential ankyrin 1 (TRPA1) antagonists.²

References

1. Revuelto, A., Ruiz-Santaquiteria, M., de Lucio, H., *et al.* Pyrrolopyrimidine vs imidazole-phenyl-thiazole scaffolds in nonpeptidic dimerization inhibitors of *Leishmania infantum* trypanothione reductase. *ACS Infect. Dis.* **5(6)**, 873-891 (2019).
2. Vallin, K.S., Sterky, K.J., Nyman, E., *et al.* N-1-Alkyl-2-oxo-2-aryl amides as novel antagonists of the TRPA1 receptor. *Bioorg. Med. Chem. Lett.* **22(17)**, 5485-5492 (2012).

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, 05/21/2020

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