

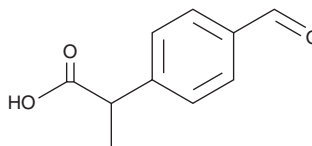
PRODUCT INFORMATION



2-(4-Formylphenyl)propionic Acid

Item No. 45718

CAS Registry No.: 43153-07-7
Formal Name: 4-formyl- α -methyl-benzeneacetic acid
Synonym: Ibuprofen Impurity K
MF: $C_{10}H_{10}O_3$
FW: 178.2
Purity: $\geq 95\%$
Supplied as: A 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

2-(4-Formylphenyl)propionic acid is supplied as a solid. A stock solution may be made by dissolving the 2-(4-formylphenyl)propionic acid in the solvent of choice, which should be purged with an inert gas. 2-(4-Formylphenyl)propionic acid is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

2-(4-Formylphenyl)propionic acid is a degradation product of ibuprofen (Item No. 70280).^{1,2}

References

1. Caviglioli, G., Valeria, P., Brunella, P., *et al.* Identification of degradation products of ibuprofen arising from oxidative and thermal treatments. *J. Pharm. Biomed. Anal.* **30(3)**, 499-509 (2002).
2. Illés, E., Takács, E., Dombi, A., *et al.* Hydroxyl radical induced degradation of ibuprofen. *Sci. Total Environ.* **447**, 286-292 (2013).

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, 05/26/2026

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