

Ibuprofen

COX inhibitor

Selective cyclooxygenase (COX) inhibitor ($IC_{50}=14.9\text{ }\mu\text{M}$). Inhibits PGH synthase-1 and PGH synthase-2 with comparable potency ($ID_{50}=8.9-14$ and $7.2-8.2\text{ }\mu\text{M}$ respectively; data are for (S) isomer). Clinically useful NSAID. Inhibits the neutrophil-dependent edema response of rabbit skin to $C5a/PGE_2$.

Citations: 4

[View Online »](#)

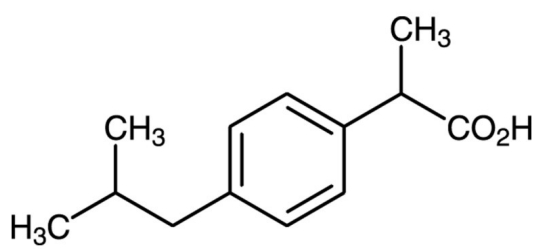
Ordering Information

[Order Online »](#)

| | |
|----------------|----|
| BML-EI164-0005 | 5g |
|----------------|----|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C. |
| Long Term Storage | Ambient |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|---|
| Alternative Name | α -Methyl-4-(2-methylpropyl)benzeneacetic acid, (\pm)-2-(4-Isobutylphenyl)propionic acid |
| Appearance | White solid. |
| CAS | 15687-27-1 |
| Couple Target | COX |
| Couple Type | Inhibitor |
| Formula | $C_{13}H_{18}O_2$ |
| MI | 14: 4881 |
| MW | 206.3 |
| Purity | $\geq 99\%$ (HPLC) |
| Solubility | Soluble in ethanol (25mg/ml). |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)