

PD 184,352

Potent and selective inhibitor of MAPK (ERK kinase 1; MEK1) activation (IC_{50} =300nM *in vitro*, IC_{50} =2nM *in vivo*). Suppresses activation of MAPK but does not block its activity. Strongly suppresses growth of human implanted colon tumors in mice.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-----|
| ALX-270-471-M001 | 1mg |
|------------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-(2-Chloro-4-iodo-phenylamino)-N-cyclopropylmethoxy-3,4-difluorobenzamide, CI-1040

Appearance White to off-white solid.

CAS 212631-79-3

Couple Target MEK

Couple Type Inhibitor

Formula $C_{17}H_{14}ClF_2IN_2O_2$

Identity Determined by NMR.

MW 478.7

Purity ≥95% (HPLC)

Solubility Soluble in organic solvents.

Source Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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