

AS-252424

PI3 kinase p110 γ inhibitor

Potent and selective phosphoinositide 3-kinase (PI(3)K) p110 γ inhibitor shown to be selective for class IB PI(3)K-mediated cellular effects.

Citations: 3

[View Online »](#)

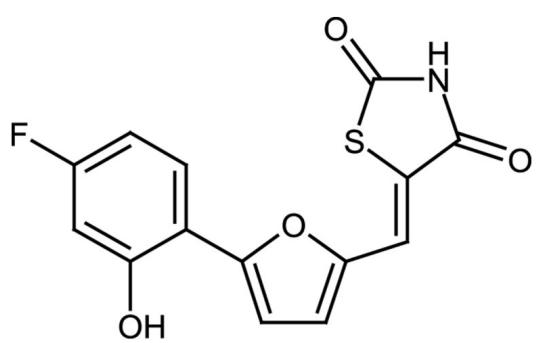
Ordering Information

[Order Online »](#)

| | |
|------------------|-----|
| ALX-270-465-M001 | 1mg |
|------------------|-----|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------|--|
| Alternative Name | 5-[5-(4-Fluoro-2-hydroxyphenyl)-furan-2-ylmethylene]-thiazolidine-2,4-dione |
| Appearance | Yellow-green to brownish solid. |
| CAS | 900515-16-4 |
| Couple Target | PI3 kinase |
| Couple Type | Inhibitor |
| Formula | $C_{14}H_8FNO_4S$ |
| MW | 305.3 |
| Purity | ≥98% (1H-NMR) |
| Solubility | Soluble in 100% ethanol (10mg/ml), DMSO (20mg/ml) or dimethyl formamide; also soluble in a 1:1 solution of DMSO:PBS, pH 7.4, (0.5mg/ml). |

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