

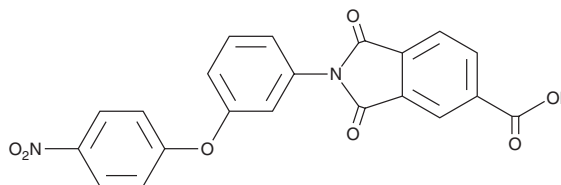
PRODUCT INFORMATION



H2L 5765834

Item No. 45940

CAS Registry No.: 420841-84-5
Formal Name: 2,3-dihydro-2-[3-(4-nitrophenoxy)phenyl]-1,3-dioxo-1H-isoindole-5-carboxylic acid
MF: C₂₁H₁₂N₂O₇
FW: 404.3
Purity: ≥98%
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

H2L 5765834 is supplied as a solid. A stock solution may be made by dissolving the H2L 5765834 in the solvent of choice, which should be purged with an inert gas. H2L 5765834 is soluble (≥10 mg/ml) in DMSO.

Description

H2L 5765834 is an antagonist of lysophosphatidic acid receptor 1 (LPA₁), LPA₃, and LPA₅ (IC₅₀s = 94, 752, and 463 nM, respectively).¹ It is selective for these receptors over LPA₂ and LPA₄. H2L 5765834 (50 μM) reverses increases in viability induced by overexpression of the gene encoding autotaxin, also known as ectonucleotide pyrophosphatase/phosphodiesterase family member 2 (ENPP2), in MIN6 pancreatic β-cells.² It decreases mechanical hypersensitivity in an NCTC clone 2472 murine fibrosarcoma model of cancer-associated pain.³

References

1. Fells, J.I., Tsukahara, R., Liu, J., *et al.* Structure-based drug design identifies novel LPA₃ antagonists. *Bioorg. Med. Chem.* **17**(21), 7457-7464 (2009).
2. Zhou, Y., Zeng, Y., Chen, Y., *et al.* Phosphodiesterase ENPP2, which is co-upregulated in obese and pregnant mice, is essential for islet beta cell compensation during obesity. *Diabetologia* **69**(4), 1000-1020 (2026).
3. Khasabova, I.A., Khasabov, S.G., Johns, M., *et al.* Exosome-associated lysophosphatidic acid signaling contributes to cancer pain. *Pain* **164**(12), 2684-2695 (2023).

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, 06/23/2026

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