

GRI977143 (LPA₂ Agonist)

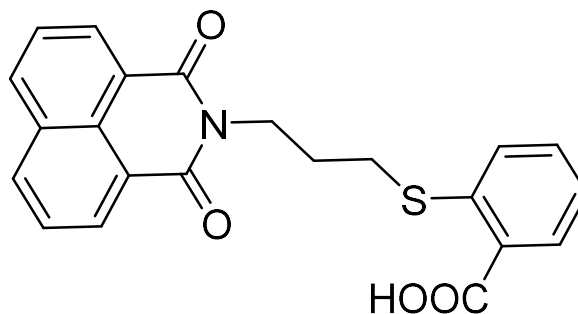
Catalog number: B-0070

Molecular Formula: C₂₂H₁₇NO₄S

MW: 391.09

CAS: 325850-81-5

Solubility: DMSO, > 5 mg/mL



Storage and Handling: Store dry at -20 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: LPA is a growth factor-like mediator and a ligand for multiple GPCR. The LPA₂ GPCR mediates antiapoptotic and mucosal barrier-protective effects in the gut. LPA GPCR activate antiapoptotic kinase pathways, inhibit apoptosis, promote cell regeneration, and augment DNA repair. GRI977143 can be used for the attenuation of programmed cell death elicited by radiation and chemotherapeutic agents, and shows strong radioprotective action by rescuing cells from apoptosis and mice irradiated with lethal doses of γ-irradiation

References: 1) Renukadevi Patil et al “Design and synthesis of sulfamoyl benzoic acid analogues with subnanomolar agonist activity specific to the LPA₂ receptor” *Journal of Medicinal chemistry*, 2014, 57, 7136–7140; 2) G.N.Kiss, et al. “Virtual screening for LPA₂-specific agonists identifies a nonlipid compound with antiapoptotic actions” *Mol. Pharmacol.* 2012, 82, 1162-1173

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.