

PRODUCT INFORMATION

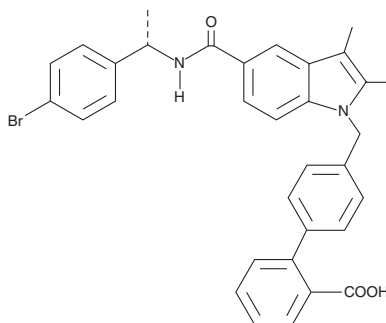


SR 1824

Item No. 11087

CAS Registry No.: 1338259-06-5
Formal Name: (S)-4'-((5-((1-(4-bromophenyl)ethyl)carbamoyl)-2,3-dimethyl-1H-indol-1-yl)methyl)-[1,1'-biphenyl]-2-carboxylic acid

MF: C₃₃H₂₉BrN₂O₃
FW: 581.5
Purity: ≥98%
UV/Vis.: λ_{max}: 252 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SR 1824 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 1824 in the solvent of choice, which should be purged with an inert gas. SR 1824 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 1824 in these solvents is approximately 30 mg/ml.

Description

Peroxisome proliferator-activated receptor γ (PPAR γ) is activated by the anti-diabetes drugs known as thiazolidinediones, including rosiglitazone (Item No. 71740) and pioglitazone (Item No. 71745).¹ Phosphorylation of PPAR γ by cyclin-dependent kinase 5 (Cdk5) causes dysregulation of genes whose expression is altered in obesity, including adiponectin.² SR 1824 is a non-agonist PPAR γ ligand ($K_i = 10$ nM) which blocks Cdk5-mediated phosphorylation.³ It does not inhibit activation of PPAR γ by rosiglitazone (Item No. 71740).³

References

1. Lehmann, J.M., Moore, L.B., Smith-Oliver, T.A., *et al.* An antidiabetic thiazolidinedione is a high affinity ligand for peroxisome proliferator-activated receptor γ (PPAR γ). *J. Biol. Chem.* **270(22)**, 12953-12956 (1995).
2. Choi, J.H., Banks, A.S., Estall, J.L., *et al.* Anti-diabetic drugs inhibit obesity-linked phosphorylation of PPAR γ by Cdk5. *Nature* **466(7305)**, 451-456 (2010).
3. Choi, J.H., Banks, A.S., Kamenecka, T.M., *et al.* Antidiabetic actions of a non-agonist PPAR γ ligand blocking Cdk5-mediated phosphorylation. *Nature* **477(7365)**, 477-481 (2011).

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, 05/22/2023

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