

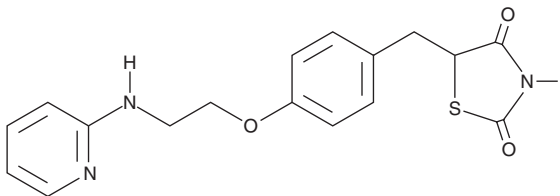
PRODUCT INFORMATION



N-desmethyl Rosiglitazone

Item No. 22103

CAS Registry No.: 257892-31-2
Formal Name: 5-[[4-[2-(2-pyridinylamino)ethoxy]phenyl]methyl]-2,4-thiazolidinedione
Synonym: SB-237216
MF: C₁₇H₁₇N₃O₃S
FW: 343.4
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 309 nm
Supplied as: A crystalline 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

N-desmethyl Rosiglitazone is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desmethyl rosiglitazone in the solvent of choice. N-desmethyl Rosiglitazone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of N-desmethyl rosiglitazone in these solvents is approximately 1, 34, and 25 mg/ml, respectively.

N-desmethyl Rosiglitazone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-desmethyl rosiglitazone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-desmethyl Rosiglitazone has a solubility of approximately 0.5 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-desmethyl Rosiglitazone is a major metabolite of rosiglitazone (Item Nos. 71740 | 71742 | 11884), a potent and selective PPAR γ ligand present in formulations that have been used to treat type 2 diabetes.^{1,2} Rosiglitazone is metabolized by the cytochrome P450 (CYP) isoform CYP2C8 to form N-desmethyl rosiglitazone.

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**, 12953-12956 (1995).
2. Kim, K.-B., Lee, D.J., Yeo, C.-W., *et al.* Simultaneous quantification of rosiglitazone and its two major metabolites, N-desmethyl and *p*-hydroxy rosiglitazone in human plasma by liquid chromatography/tandem mass spectrometry: Application to a pharmacokinetic study. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **877**(20-21), 1951-1956 (2009).

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, 11/09/2022

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