

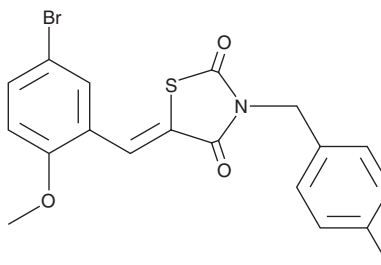
PRODUCT INFORMATION



GQ-16

Item No. 11908

CAS Registry No.: 870554-67-9
Formal Name: 5Z-[(5-bromo-2-methoxyphenyl)methylene]-3-[(4-methylphenyl)methyl]-2,4-thiazolidinedione
MF: C₁₉H₁₆BrNO₃S
FW: 418.3
Purity: ≥95%
UV/Vis.: λ_{max}: 213, 308, 359 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

GQ-16 is supplied as a crystalline solid. A stock solution may be made by dissolving the GQ-16 in the solvent of choice, which should be purged with an inert gas. GQ-16 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GQ-16 in these solvents is approximately 0.33 and 5 mg/ml, respectively.

Description

Full activation of peroxisome proliferator-activated receptor γ (PPAR γ), e.g., using thiazolidinediones, effectively treats insulin resistance and type 2 diabetes but also commonly promotes weight gain, hyperphagia, and edema.¹ GQ-16 is a partial agonist for PPAR γ (K_i = 160 nM) which induces adipogenesis significantly less than the full activator rosiglitazone (Item No. 71740).¹ Like rosiglitazone, GQ-16 reverses high fat diet-mediated impairments in insulin signaling.¹ However, GQ-16 does not induce weight gain, hyperphagia, or edema. GQ-16 does not activate PPAR α , PPAR β/δ , or RXR α .¹

Reference

1. Amato, A.A., Rajagopalan, S., Lin, J.Z., et al. GQ-16, a novel PPAR γ ligand, promotes insulin sensitization without weight gain. *J. Biol. Chem.* **287**(33), 28169-28179 (2012).

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, 03/24/2023

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