

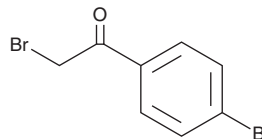
PRODUCT INFORMATION



2,4'-Dibromoacetophenone

Item No. 37358

CAS Registry No.: 99-73-0
Formal Name: 2-bromo-1-(4-bromophenyl)-ethanone
Synonyms: GSK-3 β Inhibitor VII, NSC 6224
MF: C₈H₆Br₂O
FW: 277.9
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 263 nm
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years



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

Laboratory Procedures

2,4'-Dibromoacetophenone is supplied as a solid. A stock solution may be made by dissolving the 2,4'-dibromoacetophenone in the solvent of choice, which should be purged with an inert gas. 2,4'-Dibromoacetophenone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2,4'-dibromoacetophenone in these solvents is approximately 1, 2, and 5 mg/ml, respectively.

Description

2,4'-Dibromoacetophenone is an inhibitor of glycogen synthase kinase 3 β (GSK3 β ; IC₅₀ = 0.5 μ M).¹ It is selective for GSK3 β over protein kinase A (PKA; IC₅₀ = >100 μ M). It has also been used in the derivatization of lactic acid and glycolic acid prior to analysis by HPLC.²

References

1. Conde, S., Pérez, D.I., Martínez, A., *et al.* Thienyl and phenyl α -halomethyl ketones: New inhibitors of glycogen synthase kinase (GSK-3 β) from a library of compound searching. *J. Med. Chem.* **46**(22), 4631-4633 (2003).
2. Pourasghar, M., Koenneke, A., Meiers, P., *et al.* Development of a fast and precise method for simultaneous quantification of the PLGA monomers lactic and glycolic acid by HPLC. *J. Pharm. Anal.* **9**(2), 100-107 (2019).

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, 12/21/2022

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