

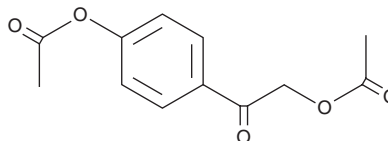
PRODUCT INFORMATION



4-(2-Acetoxyacetyl)phenyl acetate

Item No. 37888

CAS Registry No.: 42528-99-4
Formal Name: 2-(acetyloxy)-1-[4-(acetyloxy)phenyl]-ethanone
MF: $C_{12}H_{12}O_5$
FW: 236.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 250 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

4-(2-Acetoxyacetyl)phenyl acetate is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-(2-acetoxyacetyl)phenyl acetate in the solvent of choice, which should be purged with an inert gas. 4-(2-Acetoxyacetyl)phenyl acetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 4-(2-acetoxyacetyl)phenyl acetate in these solvents is approximately 1, 10, and 30 mg/ml, respectively.

Description

4-(2-Acetoxyacetyl)phenyl acetate is a synthetic intermediate that has been used in the synthesis of phenethylamine metabolites.¹

Reference

1. Holshouser, M.H. and Kolb, M. Facile synthesis of glycol metabolites of phenethylamine drugs. *J. Pharm. Sci.* **75**(6), 619-621 (1986).

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, 04/22/2025

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