PRODUCT INFORMATION



Triacetylphloroglucinol

Item No. 27073

CAS Registry No.: Formal Name:	2161-87-7 1,1',1"-(2,4,6-trihydroxy-1,3,5- benzenetriyl) <i>tris</i> -ethanone	0 0H 0
Synonym:	2,4,6-Triacetylphloroglucinol	\downarrow \downarrow \downarrow
MF:	$C_{12}H_{12}O_{6}$	
FW:	252.2	
Purity:	≥95%	но он
UV/Vis.:	λ _{max} : 273 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	ý ů
Stability:	≥4 years	

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

Laboratory Procedures

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

Description

Triacetylphloroglucinol is a C_3 -symmetric tritopic bridging ligand.¹⁻³ It has been used in the synthesis of various compounds, including trinuclear vanadium Schiff base complexes, copper complexes, and anthelmintics.

References

- 1. Spielberg, E.T. and Plass, W. 2,4,6-Triacetylphlorogucinol as threefold symmetric bridging ligand: Synthesis, structure and magnetic properties of the copper complex. Eur. J. Inorg. Chem. 20, 3093-3096 (2010).
- 2. Suresh, P., Srimurugan, S., Babu, B., et al. Asymmetric sulfoxidation of prochiral sulfides using aminoalcohol derived chiral C3-symmetric trinuclear vanadium Schiff base complexes. Tetrahedron: Asymmetry 18(23), 2820-2827 (2007).
- 3. Bowden, K., Broadbent, J.L., and Ross, W.J. Some simple anthelmintics. Br. J. Pharmacol. Chemother. 24, 714-724 (1965).

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

Suyer 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, 11/30/2022

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