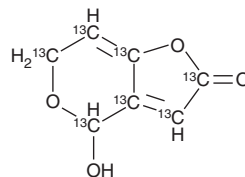


PRODUCT INFORMATION



Patulin-¹³C₇
Item No. 31301

Formal Name: 4-hydroxy-4H-furo[3,2-c]pyran-2(6H)-one-2,3,3a,4,6,7,7a-¹³C₇
Synonyms: Clairformin-¹³C₇, Expansin-¹³C₇, Mycoin C-¹³C₇
MF: [¹³C]₇H₆O₄
FW: 161.1
Purity: ≥98%
Supplied as: A 25 µg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

Patulin-¹³C₇ is intended for use as an internal standard for the quantification of patulin (Item No. 11346) by GC- or LC-MS. Patulin is a mycotoxin originally isolated from *P. griseofulvum*.¹ It is lethal to rodents (LD₅₀ = 29-55 mg/kg) and poultry (LD₅₀ = 170 mg/kg), as well as induces convulsions, dyspnea, pulmonary edema and ulceration, and gastrointestinal distension in rats, mice, chickens, and hamsters. Patulin is embryotoxic in chick eggs when used at concentrations ranging from 2.35 to 68.7 µg/egg and induces hypoplasia of the mesencephalon and telencephalon, hyperplasia of the mandibular process, and growth retardation in rat embryos. It has been found in *Penicillium*-infected apple crops.

Reference

1. Puel, O., Galtier, P., and Oswald, I.P. Biosynthesis and toxicological effects of patulin. *Toxins* **2**(4), 613-631 (2010).

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, 02/10/2026

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