# **PRODUCT** INFORMATION



Patulin-<sup>13</sup>C<sub>7</sub>

item	INO.	31	30	1

Formal Name:	4-hydroxy-4H-furo[3,2-c]pyran-2(6H)-			
	one-2,3,3a,4,6,7,7a- <sup>13</sup> C <sub>7</sub>			
Synonyms:	Clairformin- <sup>13</sup> C <sub>7</sub> , Expansin- <sup>13</sup> C <sub>7</sub> ,	H <sup>13</sup> C.		
	Mycoin $C^{-13}C_7$	H <sub>2</sub> <sup>13</sup> C 0		
MF:	[ <sup>13</sup> C] <sub>7</sub> H <sub>6</sub> O <sub>4</sub>			
FW:	161.1	0 H 13C -13C		
Purity:	≥98%	H		
Supplied as:	A 25 μg/ml solution in acetonitrile	о́н		
Storage:	-20°C			
Stability:	≥2 years			
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.				

# Description

Patulin- $^{13}C_7$  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.<sup>1</sup> It is lethal to rodents  $(LD_{50} = 29-55 \text{ mg/kg})$  and poultry  $(LD_{50} = 170 \text{ mg/kg})$ , 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.

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, 08/29/2023

# 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