PRODUCT INFORMATION



Glufosinate-d₃ (hydrochloride)

Item No. 34447

CAS Registry No.:	1323254-05-2	
Formal Name:	2-amino-4-(hydroxy(methyl-d ₃)phosphoryl)	
	butanoic acid, monohydrochloride	0
MF:	$C_5H_9D_3NO_4P \bullet HCI$	
FW:	220.6	
Chemical Purity:	≥95% (Glufosinate)	
Deuterium		
Incorporation:	≥99% deuterated forms (d ₁ -d ₃); ≤1% d ₀	• HCI
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	
Information represents	s the product specifications. Batch specific analytical resu	Ilts are provided on each certificate of analysi

Laboratory Procedures

Glufosinate-d₃ (hydrochloride) is intended for use as an internal standard for the quantification of glufosinate (Item No. 16675) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Glufosinate-d₃ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Glufosinate-d₃ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Glufosinate-d₃ (hydrochloride) is slightly soluble in methanol (heated, sonicated).

Description

Glufosinate is a broad-spectrum herbicide and racemic mixture of L-phosphinothricin, a phytotoxic compound, and D-phosphinothricin, its inactive enantiomer.¹ It is an irreversible inhibitor of glufosinate synthetase 1 (GS1) and GS2 (K.s = 1.1. and 4.8 μ M, respectively), enzymes that are required for nitrogen metabolism in plants. Glufosinate induces the production of reactive oxygen species (ROS) and lipid peroxidation in plant cell membranes. It reduces dry biomass of the weeds E. canadensis and L. rigidum with ED₅₀ values of 26 and 763 g/hectare, respectively. Glufosinate has no observed adverse effects in rats and mice when administered at doses of 4.1 and 17 mg/kg per day for three months. Formulations containing glufosinate have been used as foliar herbicides in agriculture.

Reference

1. Takano, H.K. and Dayan, F.E. Glufosinate-ammonium: A review of the current state of knowledge. Pest Manag. Sci. 76(12), 3911-3925 (2020).

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

uyer 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, 12/13/2022

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