PRODUCT INFORMATION



all-trans Retinoyl **B-D-Glucuronide**

Item No. 28057

CAS Registry No.:	401-10-5		
Formal Name:	β -D-glucopyranuronic acid, 1-retinoate		0,00
Synonyms:	Retinoyl Glucuronide,		
	Retinoyl β-D Glucuronide		,OH
MF:	$C_{26}H_{36}O_{8}$	\// o	or Y
FW:	476.6		
Purity:	≥80%		і 🔨 тон
Supplied as:	A solid		А ОН
Storage:	-80°C		on
Stability:	≥2 years		

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

Laboratory Procedures

all-trans Retinoyl β -D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the all-trans retinoyl β -D-glucuronide in the solvent of choice, which should be purged with an inert gas. all-trans Retinoyl β-D-glucuronide is slightly soluble in DMSO and methanol.

Description

all-trans Retinoyl β-D-glucuronide is a metabolite of all-trans retinoic acid (Item No. 11017) formed by the UDP-glucuronosyltransferase (UGT) system.¹ It is rapidly converted to all-trans retinoic acid following in vitro or in vivo administration.2,3

References

- 1. Salyers, K.L., Cullum, M.E., and Zile, M.H. Glucuronidation of all-trans-retinoic acid in liposomal membranes. Biochim Biophys. Acta. 11525(2), 328-334 (1993).
- 2. Rühl, R., Sass, J.O., Hau, H., et al. Effects of all-trans-retinoic acid and all-trans-retinoyl glucuronide in two in vitro systems of distinct biological complexity. Arch. Toxicol. 75(8), 497-504 (2001).
- 3. Nau, H., Elmazar, M.M.A., Rühl, R., et al. All-trans-retinoyl-β-glucuronide is a potent teratogen in the mouse because of extensive metabolism to all-trans-retinoic acid. Teratology 54(3), 150-156 (1996).

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, 07/09/2019

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