

PRODUCT INFORMATION



Recombinant Prokaryotic Lectin-Sia2 (*S. gordonii*)

Item No. 35783

Overview and Properties

Synonym:	RPL-Sia2
Source:	Recombinant His-tagged <i>S. gordonii</i> RPL-Sia2 expressed in <i>E. coli</i>
Molecular Weight:	40.72517 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥95% estimated by SDS-PAGE
Supplied in:	PBS, pH 7.2

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

Description

Recombinant prokaryotic lectin-Sia2 (RPL-Sia2) is a *S. gordonii* lectin expressed in and purified from *E. coli*.¹ It is a monomer under physiological conditions and selectively binds to glycoproteins containing terminal $\alpha(2\rightarrow3)$ -linked sialic acid (Neu5Ac) groups on O-linked glycans. RPL-Sia2 has been used to detect glycan structures in microarray experiments. RPL-Sia2 activity requires metal ions, therefore, all buffers must be supplemented with 1 mM calcium chloride, magnesium chloride, and manganese chloride. It is recommended to use metal ion-supplemented Tris-buffered saline (TBS), pH 7.6, to dilute RPL-Sia2 to the desired working concentration (≤ 20 $\mu\text{g/ml}$). Following dilution, incubate RPL-Sia2 for a minimum of 30 minutes prior to use to allow metal complexation.

Reference

1. Bertok, T., Bertokova, A., Jane, E., *et al.* Identification of whole-serum glycobiomarkers for colorectal carcinoma using reverse-phase lectin microarray. *Front. Oncol.* **11**, 735338 (2021).

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, 05/19/2022

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