

# Lectin Array 70 ab272786

### Overview

<b>Product name</b>	Lectin Array 70
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma, Cell Lysate, Tissue Lysate
<b>Assay time</b>	6h 00m
<b>Product overview</b>	Lectin Array 70 (ab272786) uses standard glass slides each spotted with 14 wells of identical lectin arrays.

This array will help researchers: 1) identify and profile the glycans in their samples; 2) determine whether their biomarker of interest has glycan moieties, and; 3) find specific glycan binding ligands in biological samples.

Lectins are glycan-binding proteins which have been purified from trees, beans and some fruits. They are highly specific for a given glycan based on their sequence and the different sugar unit structures the glycan contains. For Lectin Array 70 (ab272786) one standard glass slide is spotted with 14 identical lectin arrays, 1 in each well. Each lectin, together with the positive controls is arrayed in duplicate. The slide comes with a 16-well removable gasket which allows for the processing of 14 samples using one slide. Four slides can be nested into a tray, which matches a standard microplate and allows for automated robotic high-throughput processing of 56 arrays simultaneously. This array provides a powerful new tool for glycosylation determination, drug discovery and biomarker development; all with limited sample volumes required.

### Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	14 tests	28 tests	56 tests
20X Wash Buffer I	1 unit	1 unit	2 units
20X Wash Buffer II	1 unit	1 unit	2 units
Adhesive device sealer	2 units	4 units	8 units
Cy3 equivalent dye-conjugated Streptavidin	1 vial	2 vials	4 vials
Dialysis Vials	28 units	56 units	112 units

Components	14 tests	28 tests	56 tests
Floating Dialysis Rack	2 units	4 units	8 units
Labeling Buffer	1 vial	2 vials	4 vials
Labeling Reagent	4 vials	8 vials	16 vials
Lectin Array Glass Slide Assembly	1 slide	2 slides	4 slides
Sample Diluent	1 unit	1 unit	2 units
Slide Washer/Dryer	1 unit	1 unit	2 units
Stop Solution	1 vial	2 vials	4 vials

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors