

CUSTOMER COMMUNICATION

April 14, 2025

Re: Availability of PureLink™ Plant RNA Reagent

Dear Valued Customer,

We're writing to inform you shipment of the PureLink™ Plant RNA Reagent (Cat. No. 12322012) is delayed due to the availability of raw materials for this product. The supply issue is currently being addressed, and the estimated ship date is September 2025.

Product SKU/Code	Description
12322012	PureLink™ Plant RNA Reagent

Thank you for your continued support. Thermo Fisher Scientific is committed to supplying innovative, high-quality products for your research.

Sincerely,

Darshana Patel

Darshana Patel
Associate Product Manager, Sample Preparation