

# EUO label for instruments

Pub. No. MAN0025459 Rev. A.0

Print the following Emergency Use Only (EUO) label, then affix it on each Research Use Only (RUO) instrument authorized for use under Emergency Use Authorization (EUA). If the instrument includes labeling indicating “For Research Use Only”, cover the labeling with the EUO instrument verification label. Retain this label throughout EUA use of the instrument.

## Emergency Use Only

This instrument is authorized for use with  
Thermo Fisher Scientific assays that have  
received Emergency Use Authorization

### Revision History

Revision	Date	Description
A.0	20 May 2021	New EUO label file.

©2021 Thermo Fisher Scientific Inc. All rights reserved.

[thermofisher.com/support](https://thermofisher.com/support) | [thermofisher.com/askaquestion](https://thermofisher.com/askaquestion)  
[thermofisher.com](https://thermofisher.com)

20 May 2021

**IVD**

For In Vitro Diagnostic Use. For Emergency  
Use Authorization Only | Rx Only

**ThermoFisher**  
SCIENTIFIC