

# Stop solution 2, (50 ml)

Acid stop solution (1N HCl) intended for use as an assay stop reagent for colorimetric HRP/TMB reactions. Each bottle is sufficient for use with 5x96 wells microtiter plates. Additional assay development reagents are available as companion products to our IMMUNOSET<sup>®</sup> products.

## Ordering Information

[Order Online »](#)

|             |      |
|-------------|------|
| ADI-80-1804 | 50ml |
|-------------|------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage                      Ambient

Shipping                                      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)