

# Endonuclease ELISA Kit

A competitive immunoassay for the determination of *Serratia* endonuclease in samples

The Endonuclease (*Serratia marcescens*) ELISA is a competitive immunoassay designed to measure *Serratia* endonuclease in samples.

## Ordering Information

[Order Online »](#)

ENZ-KIT233-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Store the Standard and Detection Antibody at -20°C upon receipt. Store other kit components at 2-8° C.
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	Colorimetric detection, ELISA
Application Notes	The Endonuclease <i>Serratia marcescens</i> ELISA kit developed for quantitative determination of endonuclease in samples from downstream processing where endonuclease is used as a process or purification aid.
Assay Time	4 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Capture Secondary Antibody Microtitre coated well plate (12×8 wells) Standard (concentrated 10µg/ml) Endonuclease:HRP Conjugate Anti-Endonuclease Antibody Wash Buffer (20X) Assay Diluent TMB Substrate Stop Solution
Sensitivity	0.1 ng/ml (range 1 – 256 ng/ml)
Species Reactivity	<i>Serratia marcescens</i>
Wavelength	450 nm

Last modified: January 3, 2025



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-  
fr@enzolifesciences.com](mailto:info-fr@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)