

# Peptide YY (human) ELISA Kit

Highly sensitive Peptide YY ELISA kit  
enabling detection of PYY in serum in 3.5  
hrs.

The Peptide YY ELISA kit is a colorimetric immunometric immunoassay kit with results in 3.5 hrs.

Peptide YY (PYY) is a short 36 aa peptide released by cells in the ileum and colon in response to feeding, co-secreted with glucagon-like peptide 1 (GLP-1). In the blood and guts, PYY acts to reduce appetite. However, when injected into the CNS, PYY increases appetite. PYY plays an important role in everyday regulation of food intake.

**Manuals, SDS & CofA**

[View Online »](#)

- Sensitive measurement of PYY, detecting as little as 18.75 pg/ml
- Negligible reactivity from similar peptides
- High throughput format with results in 3.5 hrs
- Fully quantitative results that surpass semi-quantitative Western blot analysis

## Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Peptide tyrosine tyrosine, PYY, Pancreatic peptide YY
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of Peptide YY in serum, plasma, cell culture supernatant and tissue homogenate in human samples.
Assay Time	3.5 hrs
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter Plate, Standard, Sample Diluent, Detection Antibody, Wash Buffer Concentrate, Conjugate, Conjugate Diluent, Substrate, and Stop Solution
Sensitivity	18.75 pg/ml
Species Reactivity	Human
UniProt ID	P10082
Wavelength	450 nm

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)