

# 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.

## Ordering Information

[Order Online »](#)

ENZ-KIT133-0001	96 wells
-----------------	----------

## 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 [Absorbance 96 Plate Reader](#).

**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



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)