

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

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-
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)