

Total PSA (human) ELISA kit

Highly sensitive Total PSA ELISA kit enabling detection of Total PSA in human serum, plasma, urine, and tissue culture media samples in just 2.5 hours.

The Total PSA (human) ELISA kit is a colorimetric immunometric immunoassay kit with results in 2.5 hours.

Prostate-specific antigen (PSA), also known as γ -seminoprotein or kallikrein-3 (KLK3), is a kallikrein-like protease produced by the epithelial cells in the prostate gland, the lining of the urethra, and the bulbourethral gland. It is a single-chain glycoprotein with a molecular weight of about 30-33kDa in human seminal plasma and it functions normally in the liquefaction of seminal coagulum. Under healthy condition, very little PSA is secreted in the blood. PSA exists in serum in multiple forms. The proteolytic activity of PSA in blood is inhibited by irreversible formation of complexes with proteinase inhibitors such as α 1-antichymotrypsin (PSA-ACT complex), α 2-macroglobulin (not detected by immunoassays) and other acute phase proteins. A small amount of PSA is not protein bound and is called 'free PSA', which is also proteolytically inactive. Increases in glandular size and tissue damage caused by benign prostatic hypertrophy, prostatitis, or prostate cancer may increase circulating PSA levels. Monitoring the total PSA level in male blood has been used as an indicator for abnormal prostate function in the past decades. This assay provides an equal molar detection on different forms of PSA in human matrices, which improves the accuracy and sensitivity of the measurement on the total PSA levels.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

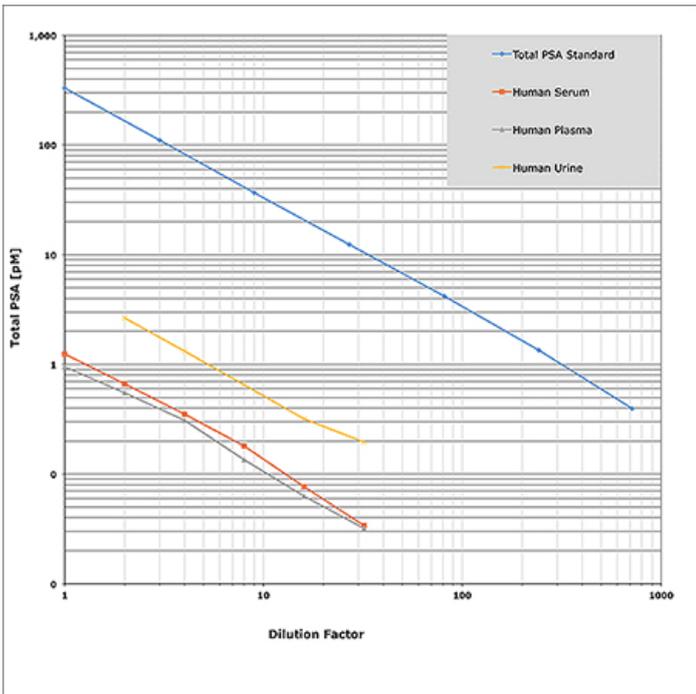
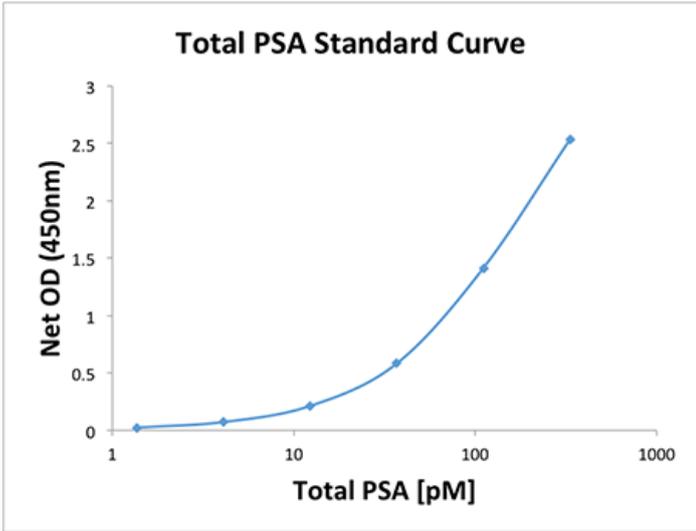
ENZ-KIT146-0001

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly-sensitive measurement of Total PSA, detecting as little as 0.287pM (or 0.024ng/ml) with negligible cross reactivity with similar proteins (KLK2, ACT)
- Compatible with various sample types (plasma, serum, urine, tissue culture media)
- High throughput format with results in 2.5 hours for up to 40 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Prostate specific antigen, Kallikrein-3
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of Total PSA in human serum, plasma, urine and tissue culture media.
Assay Time	2.5 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Microtiter plate, Assay Buffer 13, Standard, Wash Buffer Concentrate, Antibody, Antibody Diluent, Conjugate, Substrate, and Stop Solution
Crossreactivity	Minimal cross-reactivity to similar family molecules (KLK2, ACT).
Sensitivity	0.287pM or 0.024ng/ml (333-1.37pM or 28-0.12ng/ml)
Species Reactivity	Human
UniProt ID	P07288
Wavelength	450 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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