

LBP, soluble ELISA kit

The LBP, soluble ELISA kit is a complete assay for the quantitative determination of soluble LBP levels. This kit is to be used with serum, plasma and cell culture medium from human, bovine, dog, horse, porcine or rabbit origins.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-304-KI01	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	For long term storage keep vials 3 and 4 at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Lipopolysaccharide binding protein
Application	ELISA
Application Notes	For the quantitative determination of human, dog, pig, rabbit, bovine and horse soluble LBP levels in serum, plasma and cell culture medium.
Assay Time	2 hours + 10 to 12 minutes
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	1 precoated ELISA module 1 vial anti-LBP (human) detecting MAb 1 vial LBP standard (human) 1 vial reference serum 2 tablets PBS 1 vial dilution buffer 1 vial Tween 20 1 vial stopping solution 1 vial substrate solution
Quantity	96 wells (~80 tests)
Sensitivity	Range 5-50ng/ml
Species Reactivity	Bovine, Dog, Horse, Human, Porcine, Rabbit
UniProt ID	P18428 (human)
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)