



**ELISA PRODUCT INFORMATION &  
MANUAL**

**CXCL8/IL-8 ELISA Development  
Kit**

***NBP3-11747***

Enzyme-linked Immunosorbent Assay for quantitative  
detection. For research use only.

Not for diagnostic or therapeutic procedures.

[www.novusbio.com](http://www.novusbio.com) - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - [technical@novusbio.com](mailto:technical@novusbio.com)

Novus kits are guaranteed for 6 months from date of receipt

## CXCL8/IL-8 ELISA DEVELOPMENT KIT

ELISA Development Kit for quantitative determination of native or recombinant bovine CXCL8/IL-8 in solution, e.g. cell supernatant and serum/plasma samples.

<b>The kit includes</b>	<b>NBP3-11747</b> for 6 plates
Capture mAb: MT8H6 (0.5 mg/ml)	300 µl
Detection mAb: 26E5, biotinylated (0.5 mg/ml)	50 µl
Streptavidin-HRP	80 µl
Recombinant bovine IL-8 ELISA standard	1 vial
Standard reconstitution buffer A5	1 ml

To ensure total recovery of the stated quantity, vials have been overfilled.

### Shipping and storage

Shipped at ambient temperature. All reagents should be stored at 4-8 °C upon receipt, except the standard which should be stored at -20 °C. Antibodies are supplied in sterile-filtered PBS with sodium azide (0.02%). Streptavidin-HRP is supplied in PBS with 0.002% Kathon CG. The expiry date indicates how long unopened products, stored according to instructions, are recommended for use.

# General and Preparations

## Specificity

The kit contains a matched pair of monoclonal antibodies (mAbs) specific for native and recombinant human CXCL8/IL-8. The mAbs cross-react with CXCL8/IL-8 from cow, monkey and dog.

## Standard range

8-800 pg/ml

## Calibration

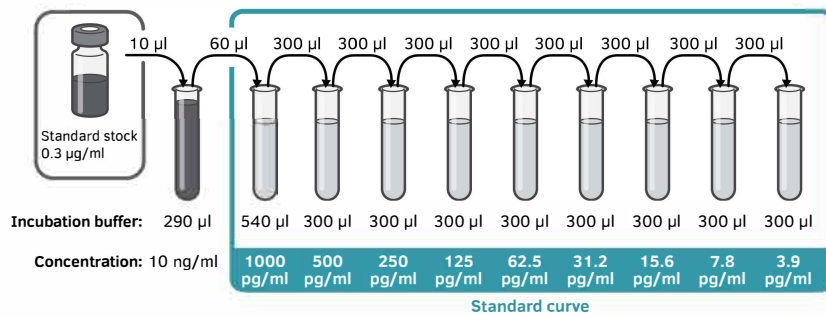
No international standard exists for calibration.

## Reconstitution of ELISA standard

Reconstitute the ELISA standard to a stock solution of 0.3 µg/ml by adding 1 ml of the standard reconstitution buffer. Allow the standard to dissolve for 5 minutes and mix thoroughly. The standard should be kept in aliquots at -20 °C. Avoid repeated freeze-thaw cycles.

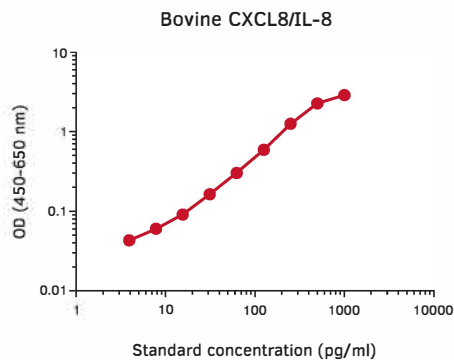
## Preparation of standard curve

Prepare within 30 minutes of use. Volumes are sufficient for duplicates.



# Protocol

- Day 1**
1. Add 100  $\mu\text{l}$ /well of capture mAb MT8H6 diluted to 2  $\mu\text{g}/\text{ml}$  in PBS, pH 7.4. Use high protein binding ELISA plates. Incubate overnight at 4-8  $^{\circ}\text{C}$ .
- Day 2**
2. Empty the plate and add 200  $\mu\text{l}$ /well of PBS with 0.05% Tween 20 and 0.1% BSA (incubation buffer) to block the plate. Incubate for 1 hour at room temperature.
  3. Wash the plate 5 times with PBS containing 0.05% Tween 20 (300  $\mu\text{l}$ /well).
  4. Add 100  $\mu\text{l}$ /well of samples or standards diluted in incubation buffer. Include assay background control, i.e. wells without standard. Incubate for 2 hours at room temperature.
  5. Wash as above.
  6. Add 100  $\mu\text{l}$ /well of detection mAb 26E5-biotin diluted to 0.1  $\mu\text{g}/\text{ml}$  in incubation buffer. Incubate for 1 hour at room temperature.
  7. Wash as above.
  8. Add 100  $\mu\text{l}$ /well of Streptavidin-HRP diluted 1:1000 in incubation buffer. Incubate for 1 hour at room temperature. Please note that sodium azide used in buffers will inhibit HRP activity.
  9. Wash as above.
  10. Add 100  $\mu\text{l}$ /well of TMB substrate and incubate for 15 minutes.
  11. Add 100  $\mu\text{l}$ /well of 0.2 M  $\text{H}_2\text{SO}_4$  to stop the reaction.
  12. Measure the optical density in an ELISA reader at 450 nm within 15 min. Preferably use a reader capable of subtracting a reference wavelength of between 570 and 650 nm. Representative standard curve shown below.



Quality management system complies with the standards  
ISO 9001:2015 & ISO 13485:2016.



The products are for research use only.

---