



ELISA PRODUCT INFORMATION & MANUAL

IL-17/IL-17A ELISA Development Kit

NBP3-11752

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

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

IL-17/IL-17A ELISA DEVELOPMENT KIT

ELISA Development Kit for quantitative determination of native goat IL-17A in solution, e.g. cell supernatant.

The kit includes		NBP3-11752 for 6 plates	
Capture mAb: MT49A7 (0.5 mg/ml)		300 µl	
Detection mAb: MT51B8, biotinylated (0.5 mg/ml)		150 µl	
Streptavidin-HRP		80 µl	
Recombinant bovine IL-17A 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 bovineIL-17/IL-17A. The mAbs cross-react with **IL-17/IL-17A** from goat. The ELISA standard is recombinant bovine IL-17/IL-17A.

Standard range

1-200 pg/ml

Calibration

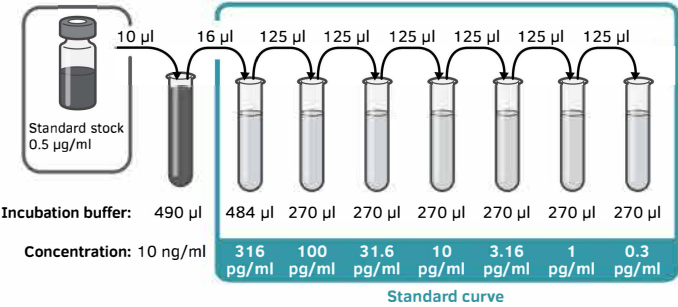
No international standard exists for calibration.

Reconstitution of ELISA standard

Reconstitute the ELISA standard to a stock solution of 0.5 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.

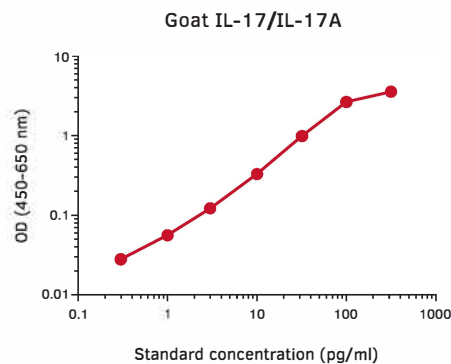
Preparation of standard curve

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



Protocol

- Day 1**
1. Add 100 µl/well of capture mAb MT49A7 diluted to 2 µg/ml in PBS, pH 7.4. Use high protein binding ELISA plates. Incubate overnight at 4-8 °C.
- Day 2**
2. Empty the plate and add 200 µl/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 µl/well).
 4. Add 100 µl/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 µl/well of detection mAb MT51B8-biotin diluted to 1 µg/ml in incubation buffer. Incubate for 1 hour at room temperature.
 7. Wash as above.
 8. Add 100 µl/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 µl/well of TMB substrate and incubate for 15 minutes.
 11. Add 100 µl/well of 0.2 M H₂SO₄ 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.

