

# Product Datasheet

## NFkB Secreted Alkaline Phosphatase Reporter Assay Kit (Colorimetric) NBP2-25286

Unit Size: 1 Kit

Storage of components varies. See protocol for specific instructions.

[www.novusbio.com](http://www.novusbio.com)



[technical@novusbio.com](mailto:technical@novusbio.com)

**Publications: 13**

Protocols, Publications, Related Products, Reviews, Research Tools and Images at:  
[www.novusbio.com/NBP2-25286](http://www.novusbio.com/NBP2-25286)

Updated 4/6/2021 v.20.1

Earn rewards for product  
reviews and publications.

Submit a publication at [www.novusbio.com/publications](http://www.novusbio.com/publications)

Submit a review at [www.novusbio.com/reviews/destination/NBP2-25286](http://www.novusbio.com/reviews/destination/NBP2-25286)



**NBP2-25286****NFκB Secreted Alkaline Phosphatase Reporter Assay Kit (Colorimetric)**

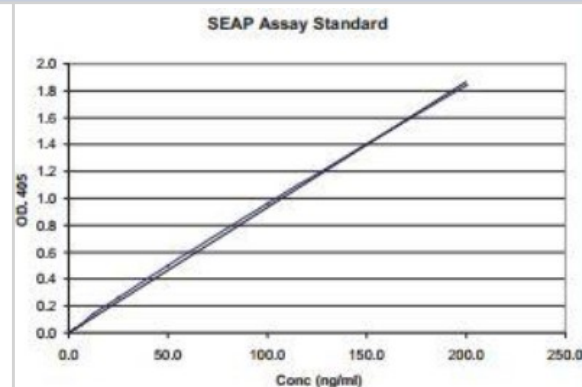
<b>Product Information</b>	
<b>Unit Size</b>	1 Kit
<b>Concentration</b>	Concentration is not relevant for this product. Please see the protocols for proper use of this product.
<b>Storage</b>	Storage of components varies. See protocol for specific instructions.

<b>Product Description</b>	
<b>Description</b>	<p>SEAP is secreted into the media so no cell lysis is required for the detection of its enzymatic activity.</p> <p>A secreted indicator function permits multiple kinetics experiments using only one culture through sequential sampling of the medium, and also allows the cells to be used for other purposes, such as RNA extraction, Western blot analysis, and other assays</p> <p>SEAP is stable in serum, allowing it to be used as a reporter assay in vivo.</p> <p>SEAP catalyzes the hydrolysis of pNitrophenyl phosphate (PNPP) producing a yellow product that can be read in a spectrophotometer or ELISA reader at 405 nm.</p>
<b>Species</b>	Human
<b>Kit Components</b>	SEAP Standard (200 ug/mL), PNPP Substrate (8 x 5 mg), pNF-kappa B/SEAP plasmid at 0.5 ug/uL (2 x 10 ug), pCMV/SEAP plasmid at 0.5 ug/uL (2 x 10 ug), 10X Sample Dilution Buffer (10 mL), 10X PNPP Buffer (4 mL), 96-well Microplate (4 EA), Microplate seal (4 EA)

<b>Product Application Details</b>	
<b>Applications</b>	Enzyme Activity
<b>Recommended Dilutions</b>	Enzyme Activity

**Images**

Enzyme Activity: NFκB Secreted Alkaline Phosphatase Reporter Assay Kit (Colorimetric) [NBP2-25286] - Standard curve for the SEAP protein provided in the NF-KB SEAPorter(TM) Assay Kit. A serial dilution of SEAP protein was added to a 96-well microtiter plate and the standard curve was generated according to the Kit assay protocol.



## Publications

Wang RW, ViganO S, Ben-David U et al. Aneuploid senescent cells activate NF-kappa B to promote their immune clearance by NK cells EMBO reports Jun 8 2021 12:00AM [PMID: 34105235]

Wang R, ViganO S, Ben-David U et al. Aneuploid cells activate NF- kappa B to promote their immune clearance by NK cells bioRxiv Jun 28 2020 12:00AM (EnzAct, Human)

Lai L, Yang G, Yao X et al. NLRC4 Mutation in flagellin-derived peptide CBLB502 ligand-binding domain reduces the inflammatory response but not radioprotective activity J. Radiat. Res. 2019 Oct 10 [PMID: 31599956] (Human)

Manzo Emiliano, Cutignano Adele, Pagano Dario et al. A new marine-derived sulfoglycolipid triggers dendritic cell activation and immune adjuvant response. Scientific Reports 2017 [PMID: 28740080]

Fereig RM, Abdelbaky HH, Kuroda Y, Nishikawa Y. Critical role of TLR2 in triggering protective immunity with cyclophilin entrapped in oligomannose-coated liposomes against Neospora caninum infection in mice Vaccine Feb 8 2019 12:00AM [PMID: 30660401] (Mouse)

Toshimitsu T, Ozaki S, Mochizuki J et al. Effects of Lactobacillus plantarum OLL2712 culture conditions on the anti-inflammatory activity for murine immune cells and obese and type 2 diabetic mice Appl. Environ. Microbiol. Jan 13 2017 12:00AM [PMID: 28087537]

O'Neill S, Humphries D, Tse G et al. Heat shock protein 90 inhibition abrogates TLR4-mediated NF-kB activity and reduces renal ischemia-reperfusion injury. Sci Rep. 2015 Aug 07 [PMID: 26248657] (ELISA)

Chen J, Ng MM, Chu JJ. Activation of TLR2 and TLR6 by Dengue NS1 Protein and Its Implications in the Immunopathogenesis of Dengue Virus Infection. PLoS Pathog. 2015 Jul 01 [PMID: 26226614] (ELISA)

Hayashi T, Yao S, Crain B et al. Induction of Tolerogenic Dendritic Cells by a PEGylated TLR7 Ligand for Treatment of Type 1 Diabetes PLoS ONE. 2015 Jun 16 [PMID: 26076454] (ELISA, Human)

McKim JM, Wilga PC, Pregonzer JF, Blakemore WR. The common food additive carrageenan is not a ligand for toll-like-receptor 4 (TLR4) in an HEK293-TLR4 reporter cell-line model Food Chem. Toxicol. 2015 Jan 29 [PMID: 25640528] (ELISA)

### Details:

SEAPorter™ assay kit used for measuring secretable alkaline phosphatase (SEAP) in experiments involving evaluation of the ability of different types of Carrageenan/CGN to bind/activate TLR4 signaling in Human TLR4/MD-2/CD14 SEAP - (SEAPorter™) Stable Reporter Cell Line [HEK293 cells stably transfected with a TLR4/MD-2/CD14/NF-kB/SEAP reporter construct - expresses full-length human TLR4, MD-2, CD14, and SEAP under the control of an NFkB response element].

Zurcher C, Sauter KS, Schweizer M. Pestiviral Erns blocks TLR-3-dependent IFN synthesis by LL37 complexed RNA. Veterinary Microbiology. 2014 Oct 13 [PMID: 25457366] (ELISA )

Signorino G, Mohammadi N, Patane F et al. Role of TLR13 in innate immune recognition of group B streptococci. Infect. Immun. 2014 Sep 15 [PMID: 25225249] (ELISA, Human)

### Details:

Human NFkB Secreted Alkaline Phosphatase SEAP - (SEAPorter™) Stable Reporter Cell Line was transiently transfected with full-length mouse TLR13 and stimulated with heat-killed bacteria and bacterial RNA or with phorbol myristate acetate /PMA followed by quantification of NFkB Secreted Alkaline Phosphatase using ELISA kit for the latter.

More publications at <http://www.novusbio.com/NBP2-25286>



## Procedures

### MSDS (NBP2-25286)

NFkB Secreted Alkaline Phosphatase Reporter Assay Kit (Colorimetric): [https://www.novusbio.com/products/nfkb-secreted-alkaline-phosphatase-reporter-kit\\_nbp2-25286](https://www.novusbio.com/products/nfkb-secreted-alkaline-phosphatase-reporter-kit_nbp2-25286)

#### Product and Contents

Product Name Nf-kB SEAPorter Assay Kit, SEAPorter Assay Kit

Catalog No. NBP2-25286

Contents p-Nitrophenyl Phosphate (KC-103), Secreted Placental Alkaline Phosphatase (IMK-515-1), SEAP Plasmids (IMK-515-3, IMK-515-4), Sodium Azide (Sample Dilution Buffer, IMK-515-6)

CAS#: 33013-2 (pNPP, KC-103)

CAS#: None (SEAP, IMK-515-1)

CAS#: None (SEAP Plasmids IMK-515-3, IMK515-4)

CAS# 26628-22-8 (10X Sample Dilution Buffer with 0.05% Sodium Azide, IMK-515-6)

p-Nitrophenyl Phosphate (pNPP) KC-103

#### Hazardous Information

Eye: Causes eye irritation.

Skin: Causes skin irritation. May be harmful if absorbed through skin.

Ingestion: May be harmful if swallowed. Call a physician

Inhalation: May causes respiratory tract irritation. May be harmful if inhaled.

Symptoms of Exposure: Irritation

#### Handling

Use with adequate ventilation. Minimize dust generation and accumulation. Avoid contact with eyes, skin, and clothing. Avoid ingestion and inhalation.

#### Storage

Store at -20 degrees C. Store in a tightly closed container. Light and moisture sensitive. Stable under recommended storage conditions

#### Personal Protection

Wear appropriate protective clothing. Protect the skin by handling with gloves. Protect eyes by wearing face shield and safety glasses. Wash hands immediately before and after use.

#### First Aid Measure

Eyes: Immediately flush eyes with plenty of water for at least 15 minutes, occasionally lifting the upper and lower eyelids. Consult a physician.

Skin: Flush skin with excess amounts of water for at least 15 minutes. Consult a physician

Ingestion: Toxic material. If swallowed, get medical aid immediately. Only induce vomiting if directed to do so by medical personnel. Never give anything by mouth to an unconscious person.

Inhalation: Remove from exposure and move to fresh air immediately. If breathing is difficult, give oxygen. **SPEED IS ESSENTIAL, OBTAIN MEDICAL AID IMMEDIATELY.** Do not use mouth-to-mouth resuscitation if victim ingested or enhaled the substance; induce artificial respiration with the aid of a pocket mask equipped with a one-way valve or other proper respiratory medical device.

Notes: Consult a physician. Show this safety data sheet to the physician. Move away from the dangerous area.

#### Secreted Alkaline Phosphatase

##### Hazardous Information

Eye: Causes eye irritation.

Skin: Causes skin irritation. May be harmful if absorbed through skin.

Ingestion: May be harmful if swallowed. Call a physician

Inhalation: May causes respiratory tract irritation. May be harmful if inhaled.

Symptoms of Exposure: Irritation

#### Handling

Handle as if capable of transmitting infectious agents. Use with adequate ventilation. Avoid contact with eyes, skin, and clothing. Avoid ingestion and inhalation.

### Storage

Store at -20 degrees C. Store in a tightly closed container. Light and moisture sensitive. Stable under recommended storage conditions

### Personal Protection

Wear appropriate protective clothing. Protect the skin by handling with gloves. Protect eyes by wearing face shield and safety glasses. Wash hands immediately before and after use.

### First Aid Measure

Eyes: Immediately flush eyes with plenty of water for at least 15 minutes, occasionally lifting the upper and lower eyelids. Consult a physician.

Skin: Flush skin with excess amounts of water for at least 15 minutes. Consult a physician

Ingestion: POISON material. If swallowed, get medical aid immediately. Only induce vomiting if directed to do so by medical personnel. Never give anything by mouth to an unconscious person.

Inhalation: Remove from exposure and move to fresh air immediately. If breathing is difficult, give oxygen. **SPEED IS ESSENTIAL, OBTAIN MEDICAL AID IMMEDIATELY.** Do not use mouth-to-mouth resuscitation if victim ingested or inhaled the substance; induce artificial respiration with the aid of a pocket mask equipped with a one-way valve or other proper respiratory medical device.

Notes: Consult a physician. Show this safety data sheet to the physician. Move away from the dangerous area.

### SEAP Plasmids

#### Hazardous Information

Eye: Causes eye irritation.

Skin: Causes skin irritation. May be harmful if absorbed through skin.

Ingestion: May be harmful if swallowed. Call a physician

Inhalation: May causes respiratory tract irritation. May be harmful if inhaled.

Symptoms of Exposure: Irritation

### Handling

Handle as if capable of transmitting infectious agents. Use with adequate ventilation. Avoid contact with eyes, skin, and clothing. Avoid ingestion and inhalation.

### Storage

Store at -20 degrees C. Store in a tightly closed container. Light and moisture sensitive. Stable under recommended storage conditions

### Personal Protection

Wear appropriate protective clothing. Protect the skin by handling with gloves. Protect eyes by wearing face shield and safety glasses. Wash hands immediately before and after use.

### First Aid Measure

Eyes: Immediately flush eyes with plenty of water for at least 15 minutes, occasionally lifting the upper and lower eyelids. Consult a physician.

Skin: Get medical aid immediately. Immediately flush skin with plenty of water for at least 15 minutes while removing contaminated clothing and shoes. **SPEEDY ACTION IS CRITICAL!**

Ingestion: POISON material. If swallowed, get medical aid immediately. Only induce vomiting if directed to do so by medical personnel. Never give anything by mouth to an unconscious person.

Inhalation: Remove from exposure and move to fresh air immediately. If breathing is difficult, give oxygen. **SPEED IS ESSENTIAL, OBTAIN MEDICAL AID IMMEDIATELY.** Do not use mouth-to-mouth resuscitation if victim ingested or inhaled the substance; induce artificial respiration with the aid of a pocket mask equipped with a one-way valve or other proper respiratory medical device.

Notes: Consult a physician. Show this safety data sheet to the physician. Move away from the dangerous area.

### 10X Sample Dilution Buffer with 0.05% Sodium Azide

#### Hazardous Information

Eye: Causes eye irritation.

Skin: Causes skin irritation. May be harmful if absorbed through skin.

Ingestion: May be fatal if swallowed.

Inhalation: May causes respiratory tract irritation. May be harmful if inhaled.

Symptoms of Exposure: Irritation

#### Handling

Use with adequate ventilation. Minimize dust generation and accumulation. Avoid contact with eyes, skin, and clothing. Avoid ingestion and inhalation.

#### Storage

Store at 4 degrees C. Store in a tightly closed container. Light and moisture sensitive. Stable under recommended storage conditions

#### Personal Protection

Wear appropriate protective clothing. Protect the skin by handling with gloves. Protect eyes by wearing face shield and safety glasses. Wash hands immediately before and after use.

#### First Aid Measure

Eyes: Immediately flush eyes with plenty of water for at least 15 minutes, occasionally lifting the upper and lower eyelids. Consult a physician.

Skin: Get medical aid immediately. Immediately flush skin with plenty of water for at least 15 minutes while removing contaminated clothing and shoes. **SPEEDY ACTION IS CRITICAL!**

Ingestion: **POISON** material. If swallowed, get medical aid immediately. Only induce vomiting if directed to do so by medical personnel. Never give anything by mouth to an unconscious person.

Inhalation: Remove from exposure and move to fresh air immediately. If breathing is difficult, give oxygen. **SPEED IS ESSENTIAL, OBTAIN MEDICAL AID IMMEDIATELY.** Do not use mouth-to-mouth resuscitation if victim ingested or enhaled the substance; induce artificial respiration with the aid of a pocket mask equipped with a one-way valve or other proper respiratory medical device.

Notes: Consult a physician. Show this safety data sheet to the physician. Move away from the dangerous area.





### **Novus Biologicals USA**

10730 E. Briarwood Avenue  
Centennial, CO 80112  
USA

Phone: 303.730.1950  
Toll Free: 1.888.506.6887  
Fax: 303.730.1966  
novus@novusbio.com

### **Bio-Techne Canada**

21 Canmotor Ave  
Toronto, ON M8Z 4E6  
Canada

Phone: 905.827.6400  
Toll Free: 855.668.8722  
Fax: 905.827.6402  
canada.inquires@bio-techne.com

### **Bio-Techne Ltd**

19 Barton Lane  
Abingdon Science Park  
Abingdon, OX14 3NB, United Kingdom

Phone: (44) (0) 1235 529449  
Free Phone: 0800 37 34 15  
Fax: (44) (0) 1235 533420  
info.EMEA@bio-techne.com

### **General Contact Information**

www.novusbio.com  
Technical Support: technical@novusbio.com  
Orders: orders@novusbio.com  
General: novus@novusbio.com

### **Limitations**

This product is for research use only and is not approved for use in humans or in clinical diagnosis. Kits are guaranteed for 6 months from date of receipt.

For more information on our 100% guarantee, please visit [www.novusbio.com/guarantee](http://www.novusbio.com/guarantee)

Earn gift cards/discounts by submitting a review: [www.novusbio.com/reviews/submit/NBP2-25286](http://www.novusbio.com/reviews/submit/NBP2-25286)

Earn gift cards/discounts by submitting a publication using this product:  
[www.novusbio.com/publications](http://www.novusbio.com/publications)

