

## SAFETY DATA SHEET

### 1. Product and Company Identification


Supplier	Novus Biologicals, LLC 10771 E Easter Ave Centennial, CO 80112, USA PHONE: 1-888-506-6887 FAX: 303-730-1966
Product Name Catalog #	Bovine IgG ELISA Kit (Colorimetric) NBP3-12512
Product Use	The Bovine IgG ELISA Kit is intended for in vitro determination of Bovine IgG in biological samples or culture media. For Research Use Only, Not for Diagnostic Use.

### 2. Hazards Identification

#### Classification of the substance or mixture

Non-toxic

#### Label Elements

Pictogram 

Signal Word: Warning

Hazard Statement: May cause skin or eye irritation

Precautionary Statement: Wash affected area with large amounts of water

Supplemental Hazard Statements: None

Other hazards: None

### 3. Composition/Information on Ingredients

The contents in the components listed as potentially hazardous in large amounts are given below:

- 1) Bovine IgG Pre-Coated 96-well Strip Plate - (<0.09% v/v) sodium azide (CAS # 26628-22-8)
- 2) Bovine IgG Standard - (<0.09% v/v) sodium azide (CAS # 26628-22-8)
- 3) Bovine IgG Detection Antibody - (0.1% v/v) Pro Clin 200™ preservative (mixture: CAS # 55965-84-9, CAS # 3251-23-8)
- 4) 20X Dilution Buffer C - no known toxicity
- 5) HRP Solution C - Buffer  
  - ≤0.013% Methylisothiazolone (CAS # 2682-20-4)
  - Streptavidin-Horseradish Peroxidase Conjugate – no known toxicity
- 6) TMB Substrate - 3,3',5,5'-tetramethylbenzidine (non-carcinogenic analog of benzidine) (CAS # 54827-17-7)
- 7) Stop Solution – 0.18 M Sulfuric Acid (CAS #7664-93-9)
- 8) 20X Wash Buffer – no known toxicity

Sodium Azide (NaN<sub>3</sub>) - Sodium azide is a hazardous chemical, however, this product is not considered to be a hazardous substance according to the OSHA Hazard Communications Standard (CFR 1910.1200), which states that, if a mixture contains less than 1% of a hazardous chemical or 0.1% of a carcinogen, the mixture shall not be considered hazardous.

Biotin - No known toxicity for this calf intestine material.

Streptavidin-Horseradish Peroxidase Conjugate - The product is not classified as dangerous according to Directive 1999/45/EC and its amendments.

Pro Clin 200™ – Liquid and mist may cause severe burns to all body tissue. May be fatal if swallowed or contacted with skin. Harmful if inhaled, affects teeth. Wash all affected areas with large volumes of water. If swallowed, give 2 glasses of water to drink and consult your physician immediately. If inhaled remove to fresh air. If necessary, give artificial respiration by mouth-to-mouth. Alternate Names: Pro-Clean 400, Kathon™ CG

Sulfuric Acid - Danger! Corrosive. Liquid and mist cause severe burns to all body tissue. May be fatal if swallowed or contacted with skin. Harmful if inhaled. Affects teeth. Cancer hazard. Strong inorganic acid mists containing sulfuric acid can cause cancer. Risk of cancer depends on duration and level of exposure.

TMB - Skin and eye irritant and suspected carcinogen.

#### 4. First Aid Measures

Oral Exposure: Wash mouth with water provided person is conscious. Do not induce vomiting. Call a physician.

Dermal Exposure: Immediately wash skin with copious amounts of water. Remove contaminated clothing and shoes. Consult a physician if irritation develops.

Eye Exposure: Flush with copious amounts of water for at least 15 minutes. Assure adequate flushing by separating the eyelids with fingers. Call a physician.

#### 5. Fire-Fighting Measures

**No specific fire or explosion hazard.**

Extinguishing Media: Water, foam, carbon dioxide, or dry chemical.

Unusual Fire & Explosion Hazard: Sodium azide forms explosive chemical compounds with lead and copper plumbing. Care must be taken to wash waste down drains with large volumes of water.

#### 6. Accidental Release Measures

If spill occurs, wear chemical resistant gloves and clean up the spill area with inert absorbent material. Avoid post-cleanup contact of chemical with skin, eyes or inhalation. Dispose of clean-up waste in suitable closed containers.

#### 7. Handling and Storage

Handling: Do not breathe vapor. Do not get in eyes, on skin, and on clothing. Avoid prolonged or repeated exposure. Wear protective clothing, chemical resistant gloves, safety goggles, and wash thoroughly after handling. No food or beverages should be carried, stored, or consumed where this material is used.

Storage: Store kit at 2 to 8 ° C. Keep tightly closed. Keep away from heat and open flame. Avoid prolonged storage at elevated temperatures as it may reduce product shelf life.

#### 8. Exposure Controls/Personal Protection

No exposure limit value known

Wash hands, forearms and face thoroughly after handling chemical products, before eating, smoking and using the lavatory and at the end of the working period.

Ventilation: No special ventilation required. Good general ventilation should be sufficient to control worker to exposure to airborne contaminants. If good general ventilation is not available then local exhaust can be used.

Protective gloves: Chemical resistant gloves

Protective eyewear: Safety eyewear that complies with an approved standard should be used in the case of liquid splashes, mists, gases, or dusts.

Other protective equipment: Lab coat to limit skin exposure. Safety shower and eyewash.

General hygiene: Wash contaminated clothing and shoes before reuse.

## 9. Physical and Chemical Properties

Physical State: Liquid, Solid

## 10. Stability and Reactivity

Stability: Product is stable under normal conditions of storage and use.

Hazardous polymerizations: Will not occur under normal conditions of storage and use.

Hazardous decomposition products: Under normal conditions and use, hazardous decomposition products should not be produced.

## 11. Toxicological Information

### United States

Acute toxicity: Not available.

Chronic toxicity: Not available.

Carcinogenicity: Not available.

Mutagenicity: Not available.

Teratogenicity: Not available.

Reproductive toxicity: Not available.

### Europe

Chronic effects: No known significant effects or critical hazards.

Carcinogenicity: No known significant effects or critical hazards.

Mutagenicity: No known significant effects or critical hazards.

Teratogenicity: No known significant effects or critical hazards.

Developmental effects: No known significant effects or critical hazards.

Fertility effects: No known significant effects or critical hazards.

## 12. Ecological Information

No known significant effects or critical hazards.

## 13. Disposal Considerations

Dispose of material in accordance with federal, state, and local regulations.

#### 14. Transportation Information

No special transport regulations

DOT:	Non-hazardous for transport
ADR (road)/RID(rail):	NA
IMDG (sea):	NA
ICAO/IATA (air):	NA

#### 15. Regulatory Information

##### U.S. Federal regulations

HCS Classification: Not regulated.

United States inventory (TSCA 8b): Not determined.

SARA 302/304/311/312 extremely hazardous substances: No products were found.

SARA 302/304 emergency planning and notification: No products were found.

SARA 302/304/311/312 hazardous chemicals: No products were found.

SARA 311/312 MSDS distribution chemical inventory – hazard identification: No products were found

Clean Water Act (CWA) 307: No products were found.

Clean Water Act (CWA) 311: No products were found.

Clean Air Act (CAA) 112 accidental release prevention: No products were found.

Clean Air Act (CAA) 112 regulated flammable substances: No products were found.

Clean Air Act (CAA) 112 regulated toxic substances: No products were found.

CEPA Toxic substances: This material is not listed.

##### Canadian Regulations

WHMIS (Canada) Not controlled under WHMIS (Canada).

Canadian ARET: This material is not listed.

Canadian NPRI: This material is not listed.

Alberta Designated Substances: This material is not listed.

Ontario Designated Substances: This material is not listed.

Quebec Designated Substances: This material is not listed.

##### EU Regulations

Risk phrases: This product is not classified according to EU legislation.

##### International Regulations:

Australia inventory (AICS): Not determined.

China inventory (IECSC): Not determined.

Korea inventory (KECI): Not determined.

Philippines inventory (PICCS): Not determined.

Japan inventory (ENCS): Not determined.

#### 16. Other Information

For research use only.

Read instructions for use before using the product. Observe the general safety regulations when handling chemicals.

The above information is believed to be correct but does not purport to be all-inclusive and is intended to be used only as a guide. Novus Biologicals, LLC shall not be liable or responsible in any way for use of either this information or the material supplied. Disposal of hazardous material may be subject to federal, state, or local laws or regulations.