

MATERIAL SAFETY DATA SHEET

SECTION1 PRODUCT AND COMPANY IDENTIFICATION

Product name:	Human TAG-72 ELISA Kit (Colorimetric) - NBP3-50054
Application	For research use only
Company:	Novus Biologicals, LLC
Address:	10771 E Easter Ave Centennial, CO 80112, USA
Email:	technical@novusbio.com
Fax:	303-730-1966
Emergency Phone:	1-888-506-6887

SECTION2 HAZARDS IDENTIFICATION

Component Items	Physical form	Hazardous Ingredient	Concentration	CAS No.
Biotinylated Detection Ab/Ag	Odorless and colorless, liquid	Proclin 300	0.04%	55965-84-9
Assay diluent	Odorless and colorless, liquid	Proclin 300	0.04%	55965-84-9
HRP Conjugate	Odorless and colorless, liquid	Proclin 300	0.04%	55965-84-9
Standard	Odorless and white/faint yellow Clear powder/ solid	Proclin 300	0.04%	55965-84-9
Substrate	Odorless and colorless, liquid	Carbamide peroxide(CP)	0.05%	124-43-6
		3,3',5,5'-tetramethylbenzidine	0.005%	54827-17-7
Stop solution	Slight pungent and colorless, liquid	Sulfuric acid (H ₂ SO ₄)	9.8%	7664-93-9

2. HAZARD STATEMENT

2.1 Proclin 300

2.1.1 Classification of the substance or mixture

Classification according to Regulation (EC) No 1272/2008 [GHS/CLP].

Sensitization, skin - Category 1

2.1.2 Label Elements

Labeling according to Regulation (EC) No 1272/2008 [GHS/CLP]

Signal Word: WARNING



Danger symbol:

Hazard Statement(s):

H317: May cause an allergic skin reaction.

Precaution Statement(s):

P261: Avoid breathing dust/fumes/gas/mist/vapours/spray.

P272: Contaminated work clothing should not be allowed out of the workplace.

P280: Wear protective gloves/protective clothing/eye protection/face protection.

P302+352: IF ON SKIN: Wash with plenty of soap and water.

P333+313: If skin irritation or rash occurs: Get medical advice/attention.

2.2 Sulfuric acid (H₂SO₄)

2.2.1 Classification of the substance or mixture

Classification according to Regulation (EC) No 1272/2008 [GHS/CLP].

Skin Corrosion/Irritation - Category 2

Eye Irritation - Category 2A

2.2.2 Label Elements

Labeling according to Regulation (EC) No 1272/2008 [GHS/CLP]

Signal Word: WARNING



Danger symbol:

Hazard Statement(s):

H315: Causes skin irritation.

H319: Causes serious eye irritation.

Precaution Statement(s):

P264: Wash hands thoroughly after handling.

P280: Wear protective gloves/protective clothing/eye protection/face protection.

P302+352: IF ON SKIN: Wash with plenty of soap and water.

P332+313: If skin irritation occurs: Get medical advice/attention.

P337+313: If eye irritation persists: Get medical advice/attention.

P305+351+338: IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

P362+364 Take off contaminated clothing and wash it before reuse.

2.3 Carbamide peroxide (CP)

2.3.1 Classification of the substance or mixture

Not a hazardous substance or mixture according to Regulation (EC) No 1272/2008.

2.3.2 Label Elements

Not a hazardous substance or mixture according to Regulation (EC) No 1272/2008.

2.4 3,3',5,5'-tetramethylbenzidine

2.4.1 Classification of the substance or mixture

Not a hazardous substance or mixture according to Regulation (EC) No 1272/2008.

2.4.2 Label Elements

Not a hazardous substance or mixture according to Regulation (EC) No 1272/2008.

SECTION3 COMPOSITION/INFORMATION ON INGREDIENTS

Ingredient	CAS No.	EC No.
Sodium chloride	7647-14-5	231-598-3
Potassium chloride	7447-40-7	231-211-8
Disodium phosphate dodecahydrate	10039-32-4	231-448-7
Potassium dihydrogen phosphate	7778-77-0	231-913-4
Tris	77-86-1	201-064-4
EDTA	60-00-4	200-449-4
Glycerol	56-81-5	200-289-5
Tween20	9005-64-5	500-018-3
Protective protein	9048-46-8	232-936-2
Mannitol	69-65-8	200-711-8
PVP40	9003-39-8	--
Proclin 300	55965-84-9	--
Carbamide peroxide(CP)	124-43-6	204-701-4
Sulfuric acid (H ₂ SO ₄)	7664-93-9	231-639-5
3,3',5,5'-tetramethylbenzidine	54827-17-7	259-364-6
N,N-Dimethylformamide (DMF)	68-12-2	200-679-5
Sodium tetraphenylborate	143-66-8	205-605-5

SECTION4 FIRST-AID MEASURES

4.1 General advice

Consult a physician. Show this safety data sheet to the doctor in attendance.

4.2 If inhaled

If breathed in, move person into fresh air. If not breathing, give artificial respiration. Consult a physician.

4.3 In case of skin contact

Take off contaminated clothing and shoes immediately. Wash off with soap and plenty of water. Consult a physician.

4.4 In case of eye contact

Rinse thoroughly with plenty of water for at least 15 minutes and consult a physician.

4.5 If swallowed

Do NOT induce vomiting. Never give anything by mouth to an unconscious person. Rinse mouth with water. Consult a physician.

SECTION5 FIRE FIGHTING MEASURES

5.1 Suitable extinguishing media

Suitable: Water spray, alcohol-resistant foam, dry chemical, carbon dioxide or appropriate foam.

For small fires, use media such as "alcohol" foam, dry chemical or carbon dioxide.

For large fires, apply water from as far as possible. Use large quantities of water applied as a mist or spray. Solid streams of water may be ineffective. Cool affected containers with flooding quantities of water.

5.2 Special precautions for fire-fighters

Protective Equipment: Wear self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

5.3 Special hazards arising from the substance or mixture

Carbon oxides, Nitrogen oxides (NOx), Sulphur oxides, Hydrogen chloride gas.

SECTION6 ACCIDENTAL RELEASE MEASURES

6.1 Person-related safety precautions

Use personal protective equipment. Avoid breathing vapors, mist or gas. Ensure adequate ventilation. Remove all sources of ignition. Evacuate personnel to safe areas. Beware of vapors accumulating to form explosive concentrations. Vapors can accumulate in low areas.

6.2 Measures for environmental protection

Prevent further leakage or spillage if safe to do so. Do not let enter drains. Discharge into the environment must be avoided.

6.3 Measures for containment and cleaning

Contain spillage, and then collect with non-combustible absorbent material (eg. sand, diatomaceous earth, vermiculite). Place in a container for disposal according to local regulations. Pick up and arrange disposal without creating dust. Sweep up and shovel. Keep in suitable, closed containers for disposal. Sweep up and shovel. Contain spillage, and then collect with an electrically protected vacuum cleaner or by wet-brushing and place in container for disposal according to local regulations (see section 13). Keep in suitable, closed containers for disposal.

SECTION7 HANDLING AND STORAGE

7.1 Handling

Wear appropriate protective clothing and safety gloves.

Avoid inhalation.

Avoid contact with eyes, skin and clothing.

Mechanical exhaust required.

Keep away from ignition sources, heat and flame.

No smoking at working site.

Incompatibilities: Strong oxidizing agents, Strong acids. Handling and unloading should be light, to prevent packaging broken, damp and cause losses.

Working place should be equipped with appropriate varieties and quantities of fire fighting equipment and leakage emergency treatment equipment.

7.2 Storage

Store in cool place. Keep container tightly closed in a dry and well-ventilated place.

Keep away from heat, sparks and flame.

Keep away from sources of ignition.

Incompatible: Strong oxidizing agents, Strong acids.

Storage place should be equipped with appropriate varieties and quantities of fire fighting equipment and leakage emergency treatment equipment.

SECTION8 EXPOSURE CONTROL/PPE

8.1 Engineering Controls

Mechanical exhaust required. Safety shower and eye bath.

8.2 Personal Protective Equipment

Respiratory: Government approved respirator if needed.

Eye/face: Chemical safety goggles if needed.

Clothing: Wear appropriate protective clothing.

Hand/skin: Protective gloves. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.

Body protection: Wear suitable protective clothing according to the concentration and amount of the substance at the workplace.

8.3 Other Protect

No smoking, drinking and eating at working site. Wash thoroughly after handling.

SECTION9 PHYSICAL/CHEMICAL PROPERTIES

9.1 Proclin 300

Appearance: Liquid

Odour: No data available

Odour threshold: No data available

pH 4.1 at 100 g/L

Melting point/freezing point: -40 °C

Initial boiling point and boiling range: 189 °C

Flash point: 118 °C - closed cup

Evaporation rate: No data available

Flammability (solid, gas): No data available

Upper/lower flammability or explosive limits: No data available

Vapour pressure: No data available

Vapour density: No data available

Relative density: 1.03 g/cm³

Water solubility: Soluble

Partition coefficient: octanol/water: No data available

Auto-ignition temperature: No data available

Decomposition temperature: No data available

Viscosity: No data available

Explosive properties: No data available

Oxidizing properties: No data available

9.2 Sulfuric acid (H₂SO₄)

Appearance: Colorless Liquid

Odor: Pungent

Odor threshold: No data available

pH: ~1

Melting point/freezing point: No data available

Boiling point/Boiling range: No data available

Flash point: No data available

Evaporation rate: No data available

Flammability (solid, gas): No data available

Upper/lower flammability or explosive limits: No data available

Vapor density: No data available

Vapor pressure: No data available

Relative density: No data available

Solubility in/Miscibility with Water: Soluble

Partition coefficient: octanol/water: No data available

Auto igniting: No data available

Decomposition temperature: No data available

Viscosity: No data available

9.3 Carbamide peroxide (CP)

Appearance: White crystalline

Odour: No data available

Odour threshold: No data available

pH: No data available

Melting point/freezing point: 90 - 93 °C - lit.

Initial boiling point and boiling range: No data available

Flash point: No data available

Evaporation rate: No data available

Flammability (solid, gas): No data available

Upper/lower flammability or explosive limits: No data available

Vapour pressure: 23.3 mmHg at 30 °C

Vapour density: No data available

Relative density: 1.390 g/cm³ at 20 °C

Water solubility: No data available

Partition coefficient: octanol/water: No data available

Auto-ignition temperature: No data available

Decomposition temperature: > 60 °C

Viscosity: No data available

Explosive properties: No data available

Oxidizing properties: The substance or mixture is classified as oxidizing with the category

Other safety information: Bulk density 0.6 - 0.7 g/L

9.4 3,3',5,5'-tetramethylbenzidine

Appearance: Liquid

Odor: No data available

Odor Threshold: No data available

pH: No data available

Melting point/freezing point: 168-171 °C - lit.

Initial boiling point and boiling range: 168 - 169 °C

Flash point: No data available
 Evaporation rate: No data available
 Flammability (solid,gas): No data available
 Upper/lower flammability or explosive limits: No data available
 Vapor pressure: No data available
 Vapor density: No data available
 Relative Density: No data available
 Water solubility: insoluble
 Partition coefficient: octanol/water: No data available
 Autoignition temperature: No data available
 Decomposition temperature: No data available
 Viscosity: No data available
 Explosive properties: No data available
 Oxidizing properties: No data available

SECTION10 STABILITY AND REACTIVITY

10.1 Reactivity

No data available

10.2 Chemical stability

Stable under recommended storage conditions

10.3 Possibility of hazardous reactions

No data available

10.4 Conditions to avoid

Heat, flames and sparks

10.5 Incompatible materials

Strong oxidizing agent, Light sensitive, Alcohols, Organic materials, Heavy metals, Powdered metals, Strong reducing agents, Amines, Mercaptans.

10.6 Hazardous decomposition products

Other decomposition products: No data available

Hazardous decomposition products formed under fire conditions: Carbon oxides, Nitrogen oxides (NOx), Sulphur oxides, Hydrogen chloride gas.

SECTION11 TOXICOLOGICAL INFORMATION

11.1 Proclin 300

Acute toxicity

LD₅₀ Oral - Rat - 862 mg/kg

LD₅₀ Dermal - Rabbit - 2,800 mg/kg

Skin corrosion/irritation

Skin - Rabbit Result: Corrosive Serious eye damage/eye irritation

Eyes - Rabbit Result: Corrosive to eyes

Respiratory or skin sensitisation - Guinea pig Result: May cause sensitisation by skin contact.

Carcinogenicity

IARC: No component of this product present at levels greater than or equal to 0.1% is identified as probable, possible or confirmed human carcinogen by IARC.

11.2 Sulfuric acid (H₂SO₄)

Acute toxicity

LD₅₀ Oral - Rat - 1530 mg/kg

LD₅₀ Dermal - Rabbit - 2730 mg/kg

LC₅₀ Inhalation- Rat - 850 mg/m³ 1 h

Skin corrosion/irritation: Can cause severe burns

Serious eye damage/irritation: Can cause severe burns

Respiratory or skin sensitization: No data available

Germ cell mutagenicity: No data available

Carcinogenicity: No data available

Reproductive toxicity: No data available

Aspiration hazard: Can cause severe burns

Ingestion: May be harmful if swallowed. Causes burns.

Skin contact: May be harmful if absorbed through skin. Causes burns.

Eye contact: Causes eye burns.

11.3 Carbamide peroxide (CP)

LD₅₀ = 4060 mg/kg (skin-rat)

Carcinogenicity:

IARC: No component of this product present at levels greater than or equal to 0.1% is identified as probable, possible or confirmed human carcinogen by IARC.

11.4 3,3',5,5'-tetramethylbenzidine

Acute toxicity

Oral: No data available

Inhalation: No data available

Dermal: No data available

Skin corrosion/irritation: No data available

Serious eye damage/eye irritation: No data available

Respiratory or skin sensitization: No data available

Germ cell mutagenicity

Test Type: Mouse

Test system: lymphocyte

Remarks: Mutation in mammalian somatic cells.

Carcinogenicity: No data available

Reproductive toxicity: No data available

Aspiration hazard: Can cause severe burns

SECTION12 ECOLOGICAL INFORMATION

12.1 Proclin 300

Ecotoxicity

No data available

Persistence and degradability

No data available

Bioaccumulative potential

No data available

Mobility in soil

No data available

Results of PBT and vPvB assessment

No data available

Other adverse effects

No data available

12.2 Sulfuric acid (H₂SO₄)

Ecotoxicity

No data available

Persistence and degradability

No data available

Bioaccumulative potential

No data available

Mobility in soil

No data available

Results of PBT and vPvB assessment

No data available

Other adverse effects

No data available

12.3 Carbamide peroxide (CP)

Ecotoxicity

No data available

Persistence and degradability

No data available

Bioaccumulative potential

No data available

Mobility in soil

No data available

Results of PBT and vPvB assessment

No data available

Other adverse effects

No data available

12.4 3,3',5,5'-tetramethylbenzidine

Ecotoxicity

No data available

Persistence and degradability

No data available

Bioaccumulative potential

No data available

Mobility in soil

No data available

Results of PBT and vPvB assessment

This substance/mixture contains no components considered to be either persistent, bioaccumulative and toxic (PBT), or very persistent and very bioaccumulative (vPvB) at levels of 0.1% or higher.

Other adverse effects

No data available

SECTION 13 DISPOSAL CONSIDERATION

13.1 Disposal methods

Dispose of waste in accordance to applicable national, regional, or local regulations. Burn in a chemical incinerator equipped with an afterburner and scrubber b highly flammable. Offer surplus and non-recyclable solutions to a licensed disposal company. Contact a licensed professional waste disposal service to dispose of this material.

13.2 Contaminated packaging

Dispose in the same manner as unused product.

SECTION 14 TRANSPORT INFORMATION

RID/ADR: Non-Hazardous for Transport: This substance is considered to be non-hazardous for transport.

IATA: Non-Hazardous for Air Transport.

IMO: Non-Hazardous for Sea Transport.

SECTION 15 REGULATORY INFORMATION

This material safety data sheet complies with the requirements of Regulation (EC) No. 1272/2008 and its amendments.

SECTION 16 OTHER INFORMATION

The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. We make no warranty of merchantability or any other warranty, express or implied, with respect to such information, and we assume no liability resulting from its use. Users should make their own investigation to determine the suitability of the information for their particular purposes. In no way shall we be liable for any claims, losses, or damages of any third party or for lost profits or any special, indirect, incidental, consequential or exemplary damages, howsoever arising from using the above information.