

Safety Data Sheet (SDS) for Enzyme-Linked Immunosorbent Assay kit

1. Identification of Substance/Mixture and of the Company

Product name: Rat COMP/Thrombospondin-5 ELISA Kit (Colorimetric)

Other means of identification: NBP2-82142

Components:

Antibody coated microwells

Standard

Detection Reagent A

Detection Reagent B

Standard Diluent (contains Proclin 300)

Assay Diluent A (contains Proclin 300)

Assay Diluent B (contains Proclin 300)

30X Wash Buffer

TMB Substrate

Stop Solution (contains Sulfuric Acid)

Recommended use: For Research Use Only.

Supplier: Novus Biologicals, LLC

Supplier Address:

10771 E Easter Ave, Centennial, CO 80112, USA

Tel: +1-888-506-6887 (Toll free) | Fax : +1-303-730-1966

Email: nb-customerservice@bio-techne.com

2. Hazard Identification:

See additional safety data sheets.

1. Identification of Substance/Mixture and of the Company

Product name: Standard Diluent (contains Proclin 300)

Recommended use: For Research Use Only.

Supplier: Novus Biologicals, LLC

Supplier Address:

10771 E Easter Ave, Centennial, CO 80112, USA

Tel: +1-888-506-6887 (Toll free) | Fax : +1-303-730-1966

Email: nb-customerservice@bio-techne.com

2. Hazard Identification:

Classification: Regulation (EC) No. 1272/2008 (CLP/EU-GHS), 29 CFR 1910.1200 (OSHA):

ProClin 300: Skin Sens. 1; H317: May cause an allergic skin reaction.

GHS Label elements, including precautionary statements

Signal Word: WARNING

Precautionary statements:

P261: Avoid breathing dust / fume / gas / mist / apours / 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.

P362+P364 Take off contaminated clothing and wash it before reuse.

Other hazards: none



3. Information on Ingredients

Description - ProClin 300

Contains	CAS No.	EC-No.	Index-No.	Content
Mixture of 5-Chloro-2-Methyl-4-Isothiazolin-3-One (26172-55-4) and 2-Methyl-4-Isothiazolin-3-One (2682-20-4)	96118-96-6	932-593-5	613-167-00-5	0.05%

4. First Aid Measures:

First aid personnel should ensure self protection.

Swallowing: Wash mouth out with water, provided person is conscious. Get medical attention.

Skin: Immediately wash skin with soap and copious amounts of water.

Inhalation: Remove to fresh air. If not breathing, perform CPR and get immediate medical attention.

Eyes: Flush the eyes with copious amounts of water for at least 15 minutes. Assure adequate flushing by separating eyelids with fingers. Consult a physician.

5. Fire Fighting Measures

Flammability: May be combustible at high temperature.

Flash points: N/A

Extinguishing media: Use dry chemical.

Special fire fighting procedures: Wear NIOSH/MSHA approved self-contained breathing apparatus and protective

clothing.

Unusual fire and explosion hazards: None known.

6. Accidental Release Measures

Person-related safety precautions: Use appropriate protective equipment to prevent contamination of skin, eyes and personal clothing.

Measures for environmental protection: Keep away from drains.

Measures for containment and cleaning: Soak up with inert absorbent material and dispose of as hazardous waste. Keep in suitable, closed containers for disposal.

7. Handling and Storage

Precautions for safe handling: Avoid contact with skin and eyes. Avoid inhalation of vapour or mist. Use normal measures for preventive fire protection.

Conditions for safe storage: Store according to product specifications.

8. Exposure Controls / Personal Protection

Control parameters: Contains no substances with occupational exposure limit values.

Individual protection measures: Wash hands thoroughly after handling chemical products and before eating, smoking or using the toilet. Appropriate techniques should be used to remove potentially contaminated clothing. Wash contaminated clothing before reusing.

Eye/face protection: Wear approved safety goggles.

Skin and body protection: Long sleeved clothing. Protective gloves..

Respiratory protection: In case of inadequate ventilation, use a suitable respirator.

9. Physical Data

Appearance: Colorless liquid

Odor: N/A

Odor threshold: N/A

pH: N/A

Melting point/range: N/A

Boiling point/Boiling range: N/A

Flash point: N/A

Evaporation rate: N/A

Flammability (solid, gas): N/A

Flammability Limits in Air: N/A

Vapor density: N/A

Vapor pressure: N/A

Solubility: N/A

Autoignition temperature: N/A

Decomposition temperature: N/A

Viscosity: N/A

10. Stability and Reactivity

Stability and Reactivity: The product is stable under recommended storage conditions.

Incompatible products: Strong acids/alkalis, strong oxidizing/reducing agents.

Conditions to avoid: Heat, moisture.

Hazardous decomposition products: Hazardous reactions will not occur under normal transport or storage conditions. Decomposition may occur on exposure to conditions or materials listed below.

11. Toxicological Information

Acute toxicity: Classified based on available data.

Chronic toxicity: Classified based on available data.

Carcinogenicity: Classified based on available data.

Sensitization: Classified based on available data.

Target Organ Effects: Classified based on available data.

12. Ecological Information

Ecotoxicity: The environmental impact of this product has not been fully investigated

13. Disposal Considerations

Disposal methods: Dispose of waste in accordance to applicable national, regional, or local regulations.

Contaminated packaging: Dispose in the same manner as unused product.

14. Transportation information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

Overland transportation (ADR/RID): Not regulated.

Transatlantic transportation (IMDG): Not dangerous according to the above specifications..

Air transportation (ICAO/IATA): Not regulated.

15. Regulatory information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

US Federal and State Regulations

Toxic Substances Control Act: Not applicable

SARA 313 Components: Not applicable

SARA 311/312 Hazards: Not applicable

CERCLA Reportable Quantity: Not applicable

California Proposition 65: Not applicable

16. Other information

This material is sold for in vitro use or research purposes only. It is not for any human or animal therapeutic or clinical diagnostic use.

Read instructions for use before using the product. Observe the general safety regulations when handling chemicals. Good laboratory practice is the best preventive measure to avoid hazards. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may be present unknown hazards and should be used with caution. Since our corporation cannot control the actual methods, volumes, or conditions of use, the company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein.

This information is prepared on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship.

1. Identification of Substance/Mixture and of the Company

Product name: Assay Diluent A (contains Proclin 300)

Recommended use: For Research Use Only.

Supplier: Novus Biologicals, LLC

Supplier Address:

10771 E Easter Ave, Centennial, CO 80112, USA

Tel: +1-888-506-6887 (Toll free) | Fax : +1-303-730-1966

Email: nb-customerservice@bio-techne.com

2. Hazard Identification:

Classification: Regulation (EC) No. 1272/2008 (CLP/EU-GHS), 29 CFR 1910.1200 (OSHA):

ProClin 300: Skin Sens. 1; H317: May cause an allergic skin reaction.

GHS Label elements, including precautionary statements

Signal Word: WARNING

Precautionary statements:

P261: Avoid breathing dust / fume / gas / mist / apours / 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.

P362+P364 Take off contaminated clothing and wash it before reuse.

Other hazards: none



3. Information on Ingredients

Description - ProClin 300

Contains	CAS No.	EC-No.	Index-No.	Content
Mixture of 5-Chloro-2-Methyl-4-Isothiazolin-3-One (26172-55-4) and 2-Methyl-4-Isothiazolin-3-One (2682-20-4)	96118-96-6	932-593-5	613-167-00-5	0.05%

4. First Aid Measures:

First aid personnel should ensure self protection.

Swallowing: Wash mouth out with water, provided person is conscious. Get medical attention.

Skin: Immediately wash skin with soap and copious amounts of water.

Inhalation: Remove to fresh air. If not breathing, perform CPR and get immediate medical attention.

Eyes: Flush the eyes with copious amounts of water for at least 15 minutes. Assure adequate flushing by separating eyelids with fingers. Consult a physician.

5. Fire Fighting Measures

Flammability: May be combustible at high temperature.

Flash points: N/A

Extinguishing media: Use dry chemical.

Special fire fighting procedures: Wear NIOSH/MSHA approved self-contained breathing apparatus and protective

clothing.

Unusual fire and explosion hazards: None known.

6. Accidental Release Measures

Person-related safety precautions: Use appropriate protective equipment to prevent contamination of skin, eyes and personal clothing.

Measures for environmental protection: Keep away from drains.

Measures for containment and cleaning: Soak up with inert absorbent material and dispose of as hazardous waste. Keep in suitable, closed containers for disposal.

7. Handling and Storage

Precautions for safe handling: Avoid contact with skin and eyes. Avoid inhalation of vapour or mist. Use normal measures for preventive fire protection.

Conditions for safe storage: Store according to product specifications.

8. Exposure Controls / Personal Protection

Control parameters: Contains no substances with occupational exposure limit values.

Individual protection measures: Wash hands thoroughly after handling chemical products and before eating, smoking or using the toilet. Appropriate techniques should be used to remove potentially contaminated clothing. Wash contaminated clothing before reusing.

Eye/face protection: Wear approved safety goggles.

Skin and body protection: Long sleeved clothing. Protective gloves..

Respiratory protection: In case of inadequate ventilation, use a suitable respirator.

9. Physical Data

Appearance: Colorless liquid

Odor: N/A

Odor threshold: N/A

pH: N/A

Melting point/range: N/A

Boiling point/Boiling range: N/A

Flash point: N/A

Evaporation rate: N/A

Flammability (solid, gas): N/A

Flammability Limits in Air: N/A

Vapor density: N/A

Vapor pressure: N/A

Solubility: N/A

Autoignition temperature: N/A

Decomposition temperature: N/A

Viscosity: N/A

10. Stability and Reactivity

Stability and Reactivity: The product is stable under recommended storage conditions.

Incompatible products: Strong acids/alkalis, strong oxidizing/reducing agents.

Conditions to avoid: Heat, moisture.

Hazardous decomposition products: Hazardous reactions will not occur under normal transport or storage conditions. Decomposition may occur on exposure to conditions or materials listed below.

11. Toxicological Information

Acute toxicity: Classified based on available data.

Chronic toxicity: Classified based on available data.

Carcinogenicity: Classified based on available data.

Sensitization: Classified based on available data.

Target Organ Effects: Classified based on available data.

12. Ecological Information

Ecotoxicity: The environmental impact of this product has not been fully investigated

13. Disposal Considerations

Disposal methods: Dispose of waste in accordance to applicable national, regional, or local regulations.

Contaminated packaging: Dispose in the same manner as unused product.

14. Transportation information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

Overland transportation (ADR/RID): Not regulated.

Transatlantic transportation (IMDG): Not dangerous according to the above specifications..

Air transportation (ICAO/IATA): Not regulated.

15. Regulatory information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

US Federal and State Regulations

Toxic Substances Control Act: Not applicable

SARA 313 Components: Not applicable

SARA 311/312 Hazards: Not applicable

CERCLA Reportable Quantity: Not applicable

California Proposition 65: Not applicable

16. Other information

This material is sold for in vitro use or research purposes only. It is not for any human or animal therapeutic or clinical diagnostic use.

Read instructions for use before using the product. Observe the general safety regulations when handling chemicals. Good laboratory practice is the best preventive measure to avoid hazards. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may be present unknown hazards and should be used with caution. Since our corporation cannot control the actual methods, volumes, or conditions of use, the company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein.

This information is prepared on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship.

1. Identification of Substance/Mixture and of the Company

Product name: Assay Diluent B (contains Proclin 300)

Recommended use: For Research Use Only.

Supplier: Novus Biologicals, LLC

Supplier Address:

10771 E Easter Ave, Centennial, CO 80112, USA

Tel: +1-888-506-6887 (Toll free) | Fax : +1-303-730-1966

Email: nb-customerservice@bio-techne.com

2. Hazard Identification:

Classification: Regulation (EC) No. 1272/2008 (CLP/EU-GHS), 29 CFR 1910.1200 (OSHA):

ProClin 300: Skin Sens. 1; H317: May cause an allergic skin reaction.

GHS Label elements, including precautionary statements

Signal Word: WARNING

Precautionary statements:

P261: Avoid breathing dust / fume / gas / mist / apours / 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.

P362+P364 Take off contaminated clothing and wash it before reuse.

Other hazards: none



3. Information on Ingredients

Description - ProClin 300

Contains	CAS No.	EC-No.	Index-No.	Content
Mixture of 5-Chloro-2-Methyl-4-Isothiazolin-3-One (26172-55-4) and 2-Methyl-4-Isothiazolin-3-One (2682-20-4)	96118-96-6	932-593-5	613-167-00-5	0.05%

4. First Aid Measures:

First aid personnel should ensure self protection.

Swallowing: Wash mouth out with water, provided person is conscious. Get medical attention.

Skin: Immediately wash skin with soap and copious amounts of water.

Inhalation: Remove to fresh air. If not breathing, perform CPR and get immediate medical attention.

Eyes: Flush the eyes with copious amounts of water for at least 15 minutes. Assure adequate flushing by separating eyelids with fingers. Consult a physician.

5. Fire Fighting Measures

Flammability: May be combustible at high temperature.

Flash points: N/A

Extinguishing media: Use dry chemical.

Special fire fighting procedures: Wear NIOSH/MSHA approved self-contained breathing apparatus and protective

clothing.

Unusual fire and explosion hazards: None known.

6. Accidental Release Measures

Person-related safety precautions: Use appropriate protective equipment to prevent contamination of skin, eyes and personal clothing.

Measures for environmental protection: Keep away from drains.

Measures for containment and cleaning: Soak up with inert absorbent material and dispose of as hazardous waste. Keep in suitable, closed containers for disposal.

7. Handling and Storage

Precautions for safe handling: Avoid contact with skin and eyes. Avoid inhalation of vapour or mist. Use normal measures for preventive fire protection.

Conditions for safe storage: Store according to product specifications.

8. Exposure Controls / Personal Protection

Control parameters: Contains no substances with occupational exposure limit values.

Individual protection measures: Wash hands thoroughly after handling chemical products and before eating, smoking or using the toilet. Appropriate techniques should be used to remove potentially contaminated clothing. Wash contaminated clothing before reusing.

Eye/face protection: Wear approved safety goggles.

Skin and body protection: Long sleeved clothing. Protective gloves..

Respiratory protection: In case of inadequate ventilation, use a suitable respirator.

9. Physical Data

Appearance: Colorless liquid

Odor: N/A

Odor threshold: N/A

pH: N/A

Melting point/range: N/A

Boiling point/Boiling range: N/A

Flash point: N/A

Evaporation rate: N/A

Flammability (solid, gas): N/A

Flammability Limits in Air: N/A

Vapor density: N/A

Vapor pressure: N/A

Solubility: N/A

Autoignition temperature: N/A

Decomposition temperature: N/A

Viscosity: N/A

10. Stability and Reactivity

Stability and Reactivity: The product is stable under recommended storage conditions.

Incompatible products: Strong acids/alkalis, strong oxidizing/reducing agents.

Conditions to avoid: Heat, moisture.

Hazardous decomposition products: Hazardous reactions will not occur under normal transport or storage conditions. Decomposition may occur on exposure to conditions or materials listed below.

11. Toxicological Information

Acute toxicity: Classified based on available data.

Chronic toxicity: Classified based on available data.

Carcinogenicity: Classified based on available data.

Sensitization: Classified based on available data.

Target Organ Effects: Classified based on available data.

12. Ecological Information

Ecotoxicity: The environmental impact of this product has not been fully investigated

13. Disposal Considerations

Disposal methods: Dispose of waste in accordance to applicable national, regional, or local regulations.

Contaminated packaging: Dispose in the same manner as unused product.

14. Transportation information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

Overland transportation (ADR/RID): Not regulated.

Transatlantic transportation (IMDG): Not dangerous according to the above specifications..

Air transportation (ICAO/IATA): Not regulated.

15. Regulatory information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

US Federal and State Regulations

Toxic Substances Control Act: Not applicable

SARA 313 Components: Not applicable

SARA 311/312 Hazards: Not applicable

CERCLA Reportable Quantity: Not applicable

California Proposition 65: Not applicable

16. Other information

This material is sold for in vitro use or research purposes only. It is not for any human or animal therapeutic or clinical diagnostic use.

Read instructions for use before using the product. Observe the general safety regulations when handling chemicals. Good laboratory practice is the best preventive measure to avoid hazards. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may be present unknown hazards and should be used with caution. Since our corporation cannot control the actual methods, volumes, or conditions of use, the company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein.

This information is prepared on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship.

1. Identification of Substance/Mixture and of the Company

Product name: Stop Solution (contains Sulfuric Acid)

Recommended use: For Research Use Only.

Supplier: Novus Biologicals, LLC

Supplier Address:

10771 E Easter Ave, Centennial, CO 80112, USA

Tel: +1-888-506-6887 (Toll free) | Fax : +1-303-730-1966

Email: nb-customerservice@bio-techne.com

2. Hazard Identification:

Classification: Regulation (EC) No. 1272/2008 (CLP/EU-GHS), 29 CFR 1910.1200 (OSHA):

Skin Corrosion/Irritation - Category 2

Serious eye damage/eye irritation - Category 2A

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

Signal Word: WARNING

Hazard Statements:

H315 Causes skin irritation.

H319 Causes serious eye irritation.

Precautionary statements:

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.

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

P321 Specific treatment (see Section 4).

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

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

P362+364 Take off contaminated clothing and wash it before reuse. Other hazards: none



3. Information on Ingredients

Contains	CAS No.	EC-No.	Index-No.	Classification	Conc.
Sulfuric Acid	7664-93-9	231-639-5	016-020-00-8	H315, H319	9.8%
Water	7732-18-5	231-791-2	-	-	90.2%

4. First Aid Measures:

First aid personnel should ensure self protection.

Swallowing: Do NOT induce vomiting. Never give anything by mouth to an unconscious person. Rinse mouth with water. Immediately consult a physician. Show this MSDS.

Skin: Immediately take off contaminated clothing or shoes. Wash with plenty of soap and water. Consult a physician.

Inhalation: Remove to fresh air. If not breathing, perform CPR and immediately consult a physician.

Eyes: Flush the eyes with copious amounts of water for at least 15 minutes. Assure adequate flushing by separating eyelids with fingers. Consult a physician immediately.

5. Fire Fighting Measures

Extinguishing media: Use dry chemical.

Special fire fighting procedures: Wear NIOSH/MSHA approved self-contained breathing apparatus and protective clothing.

Unusual fire and explosion hazards: In case of fire, toxic and corrosive gases may be formed.

6. Accidental Release Measures

Person-related safety precautions: Use appropriate protective equipment to prevent contamination of skin, eyes and personal clothing.

Measures for environmental protection: Keep away from drains.

Measures for containment and cleaning: Soak up with inert absorbent material and dispose of as hazardous waste. Keep in suitable, closed containers for disposal.

7. Handling and Storage

Precautions for safe handling: Avoid contact with skin and eyes. Avoid inhalation of vapour or mist. Use normal measures for preventive fire protection.

Conditions for safe storage: Store according to product specifications.

8. Exposure Controls / Personal Protection

Control parameters: Contains no substances with occupational exposure limit values.

Individual protection measures: Wash hands thoroughly after handling chemical products and before eating, smoking or using the toilet. Appropriate techniques should be used to remove potentially contaminated clothing. Wash contaminated clothing before reusing.

Eye/face protection: Wear approved safety goggles.

Skin and body protection: Long sleeved clothing. Protective gloves..

Respiratory protection: In case of inadequate ventilation, use a suitable respirator. If the respirator is the sole means of protection, use a full-face supplied air respirator. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

9. Physical Data

Appearance: Colorless liquid

Odor threshold: N/A

Melting point/range: N/A

Flash point: N/A

Flammability (solid, gas): N/A

Vapor density: N/A

Solubility: Soluble

Decomposition temperature: N/A

Odor: Pungent

pH: ~1

Boiling point/Boiling range: N/A

Evaporation rate: N/A

Flammability Limits in Air: N/A

Vapor pressure: N/A

Autoignition temperature: N/A

Viscosity: N/A

10. Stability and Reactivity

Stability and Reactivity: Contact with metals produces highly flammable hydrogen gas. Addition of water liberates excessive heat.

Conditions to avoid: Bases, Halides, Metals, Alkalis, Acetonitrile.

Materials to avoid: Most metals, oxidizers, reducers, bases, metal carbonates, cyanides, sulphides, carbides, oxides, metal acetylides, hydrides, halogens, organic or combustible materials, perchlorates, acetonitrile, permanganates, alcohols, picrates.

Hazardous decomposition products: Products formed under fire conditions: oxides of Sulphur, hydrogen gas.

11. Toxicological Information

Acute toxicity: Can cause severe burns upon contact while the vapors or mist are corrosive and can cause severe irritation or damage to the nose, throat and lungs. Ingestion of this product causes pain, nausea and vomiting and may be fatal if large doses are ingested.

Skin Corrosion: Can cause severe burns.

Serious Eye Damage: Can cause severe burns.

Aspiration Hazard: Can cause severe burns.

12. Ecological Information

Ecotoxicity: This product may affect the acidity (pH-factor) in water with risk of harmful effects to aquatic organisms.

13. Disposal Considerations

Disposal methods: Dispose of waste in accordance to applicable national, regional, or local regulations.

Contaminated packaging: Dispose in the same manner as unused product.

14. Transportation information

Classified according to the criteria of the UN Model Regulations as reflected in the IMDG Code, ADR, RID, DOT and IATA.

UN-Number: UN2796

UN proper shipping name: Battery fluid, acid or Sulfuric acid with not more than 51 percent acid

Transport hazard class(es): 8

Packaging group: II

Environmental hazards:

This product is not classified as environmentally hazardous according to the UN Model Regulations, nor a marine pollutant according to the IMDG Code.

Special precautions for users: No data available.

Additional Transport Information: No data available.

15. Regulatory information

This safety datasheet complies with the requirements of Regulation (EC) No. 453/2010.

US Federal and State Regulations

Toxic Substances Control Act: On TSCA Inventory

SARA 313 Components: Not applicable.

SARA 311/312 Hazards: Acute Health Hazard

CERCLA Reportable Quantity: 1,000 lbs

California Proposition 65: Not applicable.

16. Other information

This material is sold for in vitro use or research purposes only. It is not for any human or animal therapeutic or clinical diagnostic use.

Read instructions for use before using the product. Observe the general safety regulations when handling chemicals. Good laboratory practice is the best preventive measure to avoid hazards. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may be present unknown hazards and should be used with caution. Since our corporation cannot control the actual methods, volumes, or conditions of use, the company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein.

This information is prepared on our present knowledge. However, this shall not constitute a guarantee for any specific product features and shall not establish a legally valid contractual relationship.

Revision date: March.26.2024

Version: 1.2