

# NON-HAZARDOUS PRODUCT STATEMENT

### PElpro®-GMP Transfection Reagent

## A safety data sheet is not required for this product in accordance with Article 31 of Regulation (EC) No 1907/2006.

#### SECTION 1. IDENTIFICATION OF THE SUBSTANCE/MIXTURE AND OF THE COMPANY/UNDERTAKING

Trade name: PElpro®-GMP

Catalog number: 102000004, 102000008, 102000002, 102000011, 102000005, 102000007, 102000001, 102000010,

102000121, 102000131

Recommended use:

Reagent for the transfer of nucleic acids in vitro.

Uses advised against:

Not for direct injection in humans.

Company:

#### **POLYPLUS-TRANSFECTION SA**

75 rue Marguerite Perey 67400 ILLKIRCH FRANCE

Tel: +33 (0)3 90 40 61 80

Competent person responsible for the safety data sheet: support@polyplus-transfection.com

#### **SECTION 2. HAZARDS IDENTIFICATION**

#### 2.1. Classification of the substance or mixture

The mixture should not be considered as dangerous in accordance with Regulation (EC) 1907/2006.

The mixture should not be considered as hazardous in accordance with Regulation (EC) N°1272/2008.

Adverse physicochemical effects on human health and environment: no hazard.

#### 2.2. Label elements

Not Relevant.

#### **SECTION 3. COMPOSITION/INFORMATION ON INGREDIENTS**

Synthetic product in aqueous solution.

#### **SECTION 4. FIRST AID MEASURES**

In case of skin contact: wash with plenty of water and soap.

In case of contact with eyes: rinse immediately with plenty of water and seek medical advice.

In case of ingestion: do not under any circumstances induce vomiting. OBTAIN A MEDICAL EXAMINATION IMMEDIATELY.

In case of inhalation: remove casualty to fresh air and keep warm and at rest.

#### **SECTION 5. FIRE-FIGHTING MEASURES**

Suitable extinguishing media: use water spray, dry chemical, foam, carbon dioxide.

#### **SECTION 6. ACCIDENTAL RELEASE MEASURES**

Wear personal protection equipment.

Remove persons to safety.

Do not allow to enter into soil/subsoil. Do not allow to enter into surface water or drains.

Retain contaminated washing water and dispose it.

In case of entry into waterways, soil or drains, inform the responsible authorities.

Suitable material for taking up: absorbing material, sand.

Wash with plenty of water.

#### NON-HAZARDOUS PRODUCT STATEMENT

#### PElpro®-GMP Transfection Reagent

#### **SECTION 7. HANDLING AND STORAGE**

Avoid contact with skin and eyes, inhalation of vapours and mists.

Do not eat or drink while working.

Instructions as regards storage premises: adequately ventilated premises.

#### **SECTION 8. EXPOSURE CONTROLS/PERSONAL PROTECTION**

Protection of hands, eyes, skin and airways: not needed for normal use.

Anyway, operate according to good working practices.

#### **SECTION 9. PHYSICAL AND CHEMICAL PROPERTIES**

Appearance: liquid.

#### **SECTION 10. STABILITY AND REACTIVITY**

Stable under normal conditions.

#### **SECTION 11. TOXICOLOGICAL INFORMATION**

There is no toxicological data available on the mixture. Consider the individual concentration of each component to assess toxicological effects resulting from exposure to the mixture.

As far as we are currently aware, this mixture and its components have no properties potentially harmful to health.

#### **SECTION 12. ECOLOGICAL INFORMATION**

Adopt good working practices, so that the product is not released into the environment.

Since no ecotoxicological data about the mixture is available, the concentration of each substance must be considered to assess the ecotoxicological effects resulting from exposure to the mixture.

The product is considered to present no hazard to the environment, particularly in respect of its mobility in the soil, its persistence, its potential for bioaccumulation, its aquatic toxicity and, generally, its ecotoxicity.

#### **SECTION 13. DISPOSAL CONSIDERATIONS**

Recover if possible. In so doing, comply with the local and national regulations currently in force.