

## PEIpro<sup>®</sup>-GMP Transfection Reagent

according to Federal Register / Vol. 77, No. 58 / Monday, March 26, 2012 / Rules and Regulations

### SECTION 1. IDENTIFICATION

#### 1.1. Identification

Product form	: Mixture
Product name	: PEIpro <sup>®</sup> -GMP
Product code	: 102000004, 102000008, 102000002, 102000011, 102000005, 102000007, 102000001, 102000010, 102000121, 102000131

#### 1.2. Recommended use and restrictions on use

Use of the substance/mixture	: Reagent for the transfer of nucleic acids <i>in vitro</i> . Not for direct injection in humans.
------------------------------	--

#### 1.3. Supplier

**POLYPLUS-TRANSFECTION SA**  
75 rue Marguerite Perey  
67400 ILLKIRCH  
FRANCE  
Tel: +33 (0)3 90 40 61 80  
[support@polyplus-transfection.com](mailto:support@polyplus-transfection.com)

#### 1.4. Emergency telephone number

No additional information available

### SECTION 2. HAZARD(S) IDENTIFICATION

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

##### GHS-US classification

The mixture should not be considered as dangerous.  
The mixture should not be considered as hazardous.  
Adverse physicochemical effects on human health and environment: no hazard.

#### 2.2. GHS Label elements, including precautionary statements

##### GHS-US labelling

Hazard pictograms (GHS-US)	: No pictogram
Signal word (GHS-US)	: No signal word
Hazard statements (GHS-US)	: No statement
Precautionary statements (GHS-US)	: No statement

#### 2.3. Other hazards which do not result in classification

Other hazards not contributing to the classification	: None under normal conditions.
--	---------------------------------

#### 2.4. Unknown acute toxicity (GHS US)

Not applicable

### SECTION 3. COMPOSITION/INFORMATION ON INGREDIENTS

#### 3.1. Substances

Not applicable

#### 3.2. Mixtures

Synthetic product in aqueous solution.

## PEIpro®-GMP Transfection Reagent

## SECTION 4. FIRST AID MEASURES

## 4.1. Description of first aid measures

- First-aid measures after inhalation : Move the affected person away from the contaminated area and into the fresh air. Get medical advice/attention if you feel unwell.
- First-aid measures after skin contact : Wash with soapy water. If skin irritation or rash occurs: Get medical advice/attention.
- First-aid measures after eye contact : Immediately rinse with water for a prolonged period while holding the eyelids wide open. If eye irritation persists: Get medical advice/attention.
- First-aid measures after ingestion : Never attempt to induce vomiting. Call a physician immediately.

## 4.2. Most important symptoms and effects (acute and delayed)

- Symptoms/effects : No specific effects and/or symptoms are known.

## 4.3. Immediate medical attention and special treatment, if necessary

No immediate medical care or specific treatment identified.

## SECTION 5. FIRE-FIGHTING MEASURES

## 5.1. Suitable (and unsuitable) extinguishing media

- Suitable extinguishing media : Water spray. Water spray with additive . Carbon dioxide (CO<sub>2</sub>). Alcohol-resistant foam. Dry powder.

## 5.2. Specific hazards arising from the chemical

- Fire hazard : Flammable liquid and vapour. The vapours are denser than air and may travel along the ground. Distance ignition possible.
- Hazardous combustion products : Toxic fumes may be released Carbon oxides (CO, CO<sub>2</sub>)

## 5.3. Special protective equipment and precautions for fire-fighters

- Firefighting instructions : Cool down the containers exposed to heat with a water spray. Contain the extinguishing fluids by bunding.
- Protection during firefighting : Do not attempt to take action without suitable protective equipment. Self-contained breathing apparatus. Complete protective clothing.

## SECTION 6. ACCIDENTAL RELEASE MEASURES

## 6.1. Personal precautions, protective equipment and emergency procedures

- General measures : Only qualified personnel equipped with suitable protective equipment may intervene.

## 6.1.1. For non-emergency personnel

- Emergency procedures : Avoid contact with skin and eyes.

## 6.1.2. For emergency responders

- Protective equipment : Do not attempt to take action without suitable protective equipment.

## 6.2. Environmental precautions

Contain the spilled material by bunding. Do not allow product to spread into the environment. Do not discharge into drains or rivers. Notify authorities if product enters sewers or public waters.

## 6.3. Methods and material for containment and cleaning up

- For containment : Liquid spill: take up in sand, earth, vermiculite.

**PEIpro®-GMP Transfection Reagent**

Methods for cleaning up	: Wash non-recoverable remainder with large amounts of water.
Other information	: Dispose of contaminated materials in accordance with current regulations.

**6.4. Reference to other sections**

For further information refer to section 13.

**SECTION 7. HANDLING AND STORAGE****7.1. Precautions for safe handling**

Precautions for safe handling	: Ensure good ventilation of the work station. Earth the equipment. No flames, no sparks. Eliminate all sources of ignition. Avoid contact with skin and eyes. Do not breathe vapours. Smoking is forbidden.
Hygiene measures	: Do not drink, eat or smoke in the workplace. Wash hands and other exposed areas with mild soap and water before eating, drinking or smoking and when leaving work.

**7.2. Conditions for safe storage, including any incompatibilities**

Storage conditions	: Keep container tightly closed. Store in dry, cool, well-ventilated area. Keep away from heat. Keep away from sources of ignition. Protect from sunlight. Keep away from food, drink and animal feeding stuffs.
Incompatible products	: Combustible materials.
Packaging materials	: Original packaging.

**SECTION 8. EXPOSURE CONTROLS/PERSONAL PROTECTION****8.1. Control parameters**

Not applicable.

**8.2. Appropriate engineering controls**

Appropriate engineering controls	: Ensure good ventilation of the work station.
----------------------------------	--

**8.3. Individual protection measures/Personal protective equipment**

**Hand protection:** Not required for normal conditions of use

**Eye protection:** Not required for normal conditions of use

**Skin and body protection:** Not required for normal conditions of use

**Respiratory protection:** Not required for normal conditions of use

**SECTION 9. PHYSICAL AND CHEMICAL PROPERTIES****9.1. Information on basic physical and chemical properties**

Physical state	: Liquid
Colour	: No data available
Odour	: No data available
Odour threshold	: No data available
pH	: No data available

**PEIpro®-GMP Transfection Reagent**

Melting point	: No data available
Freezing point	: No data available
Boiling point	: No data available
Flash point	: No data available
Relative evaporation rate (butylacetate=1)	: No data available
Flammability (solid, gas)	: Not applicable.
Vapour pressure	: No data available
Relative vapour density at 20 °C	: No data available
Relative density	: No data available
Solubility	: Water: Soluble
Log Pow	: No data available
Auto-ignition temperature	: No data available
Decomposition temperature	: No data available
Viscosity, kinematic	: No data available
Viscosity, dynamic	: No data available
Explosive limits	: No data available
Explosive properties	: No data available
Oxidising properties	: No data available

**9.2. Other information**

No additional information available

**SECTION 10. STABILITY AND REACTIVITY****10.1. Reactivity**

No additional information available.

**10.2. Chemical stability**

Stable in use and storage conditions as recommended in item 7.

**10.3. Possibility of hazardous reactions**

No additional information available.

**10.4. Conditions to avoid**

Sparks. Heat. Open flame. Ignition sources.

**10.5. Incompatible materials**

Combustible materials. Alkali metals. Alkaline earth metals. Strong oxidizing agents. Strong reducing agents. Mineral acids. Peroxides. Nitrogen compounds.

**PEIpro®-GMP Transfection Reagent****10.6. Hazardous decomposition products**

Under normal conditions of storage and use, hazardous decomposition products should not be produced.

**SECTION 11. TOXICOLOGICAL INFORMATION****11.1. Information on toxicological effects**

Acute toxicity	: Not classified
Skin corrosion/irritation	: Not classified
Serious eye damage/irritation	: Not classified
Respiratory or skin sensitisation	: Not classified
Germ cell mutagenicity	: Not classified
Carcinogenicity	: Not classified
Reproductive toxicity	: Not classified
Specific target organ toxicity (single exposure)	: Not classified
Specific target organ toxicity (repeated exposure)	: Not classified
Aspiration hazard	: Not classified
Symptoms/effects	: No specific effects and/or symptoms are known.

**SECTION 12. ECOLOGICAL INFORMATION****12.1. Toxicity**

Ecology - general	: Presents no particular risk to the environment, provided the disposal requirements (see section 13) and national or local regulations are complied with.
-------------------	--

**12.2. Persistence and degradability**

Not classified

**12.3. Bioaccumulative potential**

Not classified

**12.4. Mobility in soil**

Not classified

**12.5. Other adverse effects**

Effect on global warming	: No known effects from this product.
GWPmix comment	: No known effects from this product.

**SECTION 13. DISPOSAL CONSIDERATIONS****13.1. Disposal methods**

Waste treatment methods	: Dispose of in accordance with relevant local regulations.
Additional information	: Empty the packaging completely prior to disposal. Recycle or dispose of in compliance with current legislation.

**SECTION 14. TRANSPORT INFORMATION**

Not classified

**SECTION 15. OTHER INFORMATION**

No other information