

FectoVIR[®]-AAV 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: **FectoVIR[®]-AAV GMP**
Catalog number: 102000006, 102000009, 102000012

Recommended use:
Reagent for the transfer of nucleic acids *in vitro*.
Uses advised against:
Not intended for human diagnostic or therapeutic use.
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
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.

FectoVIR®-AAV 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 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.