

pPLUS[®] AAV-Helper - Plasmid Helper

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

1.1. Product Identifier

Mixture identification: Plasmid Helper
Trade name: pPLUS[®] AAV-Helper
Catalog number: 101000181, 101000182, 101000183.

1.2. Relevant identified uses of the substance/mixture and uses advised against

Recommended use:
Plasmid Helper used for research purpose and manufacturing of viral vector.
Uses advised against:
Not for use in human.

1.3. Details of the supplier of the safety data sheet

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

1.4. Emergency telephone number

National Poisons Information Service: + 44 (0)844 892 0111
On-line information database TOXBASE (www.toxbase.org).

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 applicable.

SECTION 3. COMPOSITION/INFORMATION ON INGREDIENTS

Plasmid Helper in Tris-EDTA buffer

SECTION 4. FIRST AID MEASURES

4.1. Description of 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

pPLUS® AAV Helper - Plasmid Helper

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.

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.

End of the Non-Hazardous Product Statement.