



Safety Data Sheet

Section 1 - Product and Company Information

Product Identifier:	PLTGold®
Product Code:	PLTGold27R, PLTGold100R, PLTGold500R, PLTGold27GMP, PLTGold100GMP, PLTGold500GMP
Company Name:	Mill Creek Life Sciences, LLC
Street Address:	221 1 st Avenue SW, Suite 209, Rochester, Minnesota 55902 USA
Technical Phone:	507-287-6257
Emergency Phone:	507-287-6257
Description:	Human Platelet Lysate
Recommended Use:	Cell Culture Supplement

Section 2 - Hazard Identification

Potential Hazards:	This material is manufactured from biological materials (platelets) obtained from human donors. All apheresis platelets used in the manufacture of PLTGold® met donor eligibility criteria at the time of their collection as per applicable standards.
---------------------------	--

Section 3 - Composition/Information on Ingredients

Human platelet lysate is manufactured from donor platelets.

Section 4 - First Aid Measures

Eye Contact:	Flush immediately with water and seek medical advice.
Skin Contact:	Wash area immediately with soap and water. Remove any contaminated clothing/shoes.
Inhalation:	Remove to fresh air. If not breathing, give artificial respiration. If breathing is difficult, give oxygen. If not breathing/breathing is difficult seek medical attention.
Ingestion:	Wash mouth with water if person is conscious and seek medical advice.
Most important symptoms/ effects (acute and delayed):	See Section 11.
Indication immediate medical treatment is needed:	No data available



Section 5 - Fire Fighting Measures

Suitable extinguishing media:	Standard extinguishing media/equipment is sufficient.
Specific hazards arising from the chemical:	Use of this product in accordance with Standard Precautions should not result in flammable or explosive mixtures being formed.
Special protective equipment and precautions for fire fighters:	Standard protective equipment is sufficient.

Section 6 - Accidental Release Measures

Personal precautions, protective equipment, and emergency procedures:	See sections 7 and 8.
Methods and materials for containment and clean up:	Absorb material using common absorbent materials. Clean spill area with disinfectant used for biologic materials. Deposit waste in appropriate waste containers for removal and disposal as per Section 13.

Section 7 - Handling and Storage

Precautions for safe handling:	Follow Standard Precaution procedures.
Conditions for safe storage:	Store at -20°C or colder until ready for use. Product may be stored short term at 4°C. Keep closed when not in use.

Section 8 - Exposure Controls/Personal Protective Equipment

Exposure Limits:	Not Applicable; no exposure limit.
Appropriate engineering controls:	Ventilation is recommended for normal product usage. Area ventilation is usually adequate. If ventilation is insufficient wear suitable respiratory equipment.
Individual protection measures:	Protective gowning materials should be worn to limit exposure to eyes, skin and mucous membranes including splash resistant safety goggles/glasses, chemical resistant gloves, and impervious clothing appropriate to the quantity and usage. Dispose of contaminated gowning materials in biohazardous waste. Follow Standard Precaution procedures when using this product. Wash hands thoroughly with soap and water after using this product.

Section 9 - Physical and Chemical Properties

Appearance:	Physical state is translucent amber liquid.
Odor:	No Data Available
Odor threshold:	No Data Available



pH:	6.8-7.8
Melting/freezing point:	No Data Available
Initial boiling point/range:	No Data Available
Flash point:	No Data Available
Evaporation rate:	No Data Available
Flammability (solid, gas):	No Data Available
Upper/Lower flammability or explosive limits:	No Data Available
Vapor pressure/density:	No Data Available
Relative density:	1.021 g/mL
Solubility(ies):	Hydrophilic
Partition coefficient:	No Data Available
Auto-ignition temp.:	No Data Available
Decomposition temp.:	No Data Available
Viscosity:	Variable; gel formation may occur due to presence of clotting factors.

Section 10 - Stability and Reactivity

Reactivity:	Product is nonreactive under normal conditions of storage and use.
Stability:	Product is stable under normal conditions of storage and use.
Possibility of hazardous conditions:	Use of these products in accordance with Good Laboratory Practices should not result in hazardous reactions.
Conditions to avoid:	No Data Available
Incompatible materials:	No Data Available
Hazardous decomposition products:	No Data Available

Section 11 - Toxicological Information

Toxicological effects:	Eye Contact: No Data Available Skin Contact: No Data Available Inhalation: No Data Available Ingestion: No Data Available
Symptoms related to physical, chemical and toxicological characteristics:	No Data Available



Delayed, immediate or chronic effects of short-term/long-term exposure: No Data Available

Numeric measures of toxicity: No Data Available

Possible carcinogenic effects: No Data Available

Section 12 - Ecological Information

Ecotoxicity: No Data Available

Persistence and degradability: No Data Available

Bioaccumulative potential: No Data Available

Mobility in soil: No Data Available

Other adverse effects: No Data Available

Section 13 - Disposal Considerations

Dispose of waste and any contaminated materials in accordance with all federal, state and local environmental regulations.

Section 14 - Transport Information

This substance is considered non-hazardous for ground or air transport.

Section 15 - Regulatory Information

Not Applicable

Section 16 - Other Information

Prepared by: Mill Creek Life Sciences Quality Department

Revision Date: 06 Sep 2018

Disclaimer: The above information is believed to be correct to the best of our knowledge; it does not claim to be all inclusive and should be used as a guide only. Mill Creek shall have no liability for any direct, indirect, consequential or incidental damages arising out of the use, the results of use, the inability to use, the handling or contact with these products.

This product is intended to be used for cell culture only. Product is a reagent (not for systemic or topical use).