

# Material Safety Data Sheet

Revision Date: April 21, 2020

1. IDENTIFICATION OF	THE SUBSTRANCE/PREPARATION AND THE
	COMPANY/UNDERTAKING

Product code QP035, QP036, QP037, QP038, QP039, QP045, QP046, QP047, QP048, QP049

**Product name** BlazeTaq<sup>™</sup> Probe qPCR mix

Contact manufacturer

#### GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101 Rockville, MD 20850 USA

Phone: 301-762-0888

Fax: 301-762-3888

## 2. COMPOSITION/INFORMATION ON INGREDIENTS

#### Hazardous/Non-hazardous Components

The product contains no substances which at their given concentration, are considered to be hazardous to health. We recommend handling al chemicals with caution.

## 3. HAZARDS IDENTIFICATION

**Emergency overview** 

not applicable

Information pertaining to particular dangers for man and environment

not applicable

#### **Classification system**

The classification was made according to OSHA 29CFR 1910.1200 and 1910.1030, and is expanded upon from company and/or literature information.

4. FIRST AID MEASURES			
General information	No special measures required.		
After inhalation	Supply fresh air; consult doctor in case of complaints.		
After skin contact	Rinse with plenty of water. Generally, the product does not irritate the skin.		
After eye contact	Rinse opened eye for several minutes under running water.		
After swallowing	Not expected to present a significant ingestion hazard under anticipated conditions of normal use.		

# 5. FIRE-FIGHTING MEASURES

#### Suitable extinguishing agents

CO2, extinguishing powder or water spray. Fight larger fires with water spray or alcohol resistant foam.

Protective equipment No special measures required.

6. ACCIDENTAL REL	EASE MEASURES			
Person-related safety precautions Not required				
Measures for environmental protection Dilute with plenty of water.				
<b>Measures for cleaning/collecting</b> Absorb with liquid-binding material (sand, diatomite, acid binders, universal binders, sawdust).				
Additional information No dangerous substances are released.				
7. HANDLING AND STORAGE				
Handling Information for safe handling				
No special measures required.				
No special precautions are necessary if used correctly.				
Information about protection against explosions and fires	No special measures required.			
Storage				
Requirements to be met by storeroomsand receptacles	According to productspecification			
Information about storage in one commonstorage facility	Not required.			
Further information about storage conditions None.				

8. EXPOSURE CONTROLS / PERSONAL PROTECTION				
Additional information about design of technical systems No further data; see item 7.				
Components with limit values that require monitoring at the workplace Not required.				
Additional information	<b>n</b> The lists that were valid during the creation were used as basis.			
Personal protective equipment General protective and hygienic measures The usual precautionary measures for handling chemicals should be followed.				
Breathing equipment	Not required.			
Protection of hands	Protective gloves.			
Material of gloves	Synthetic gloves			
<b>Penetration time of glove material</b> The exact break through time has to be found out by the manufacturer of the protective gloves and has to be observed.				
Eye protection	Safety glasses			

General informationFormliquidColorColorlessOdorNo data available

Change in condition<br/>Melting point/melting range<br/>Boiling point/boiling range°C Mixture has not been tested<br/>°C Mixture has not been testedFlash point°C Mixture has not been testedDanger of explosionDo not present an explosion hazard.Vapor pressure at 20°CMixture has not been testedDensity at 20°CMixture has not been testedSolubility in / Miscibilitywith WaterFully misciblePH value at 20°C% 6

8.6
25%
75%

# 10. STABILITY AND REACTIVITY

**Thermal decomposition / conditions to be avoided** No decomposition if used according to specifications.

 Materials to be avoided

 Dangerous reactions
 No dangerous reactions known

 Dangerous products of decomposition
 No dangerous decomposition products known

# 11. TOXICOLOGICAL INFORMATION

Acute ToxicityPrimary irritant effectOn the skinNo irritant effect.On theeyeNo irritant effect.

Sensitization No sensitizing effects known.

Target organs Not applicable.

## 12. ECOLOGICAL INFORMATION

Generally not hazardous for water.

## 13. DISPOSAL CONSIDERATIONS

Dispose of in accordance with local regulations

14. TRANSPORT INFORMATION		
IATA Proper shi Hazard Cla Subsidiary Packing gi UN-No	Class	Not classified as dangerous in the meaning of transport regulations No information available No information available No information available UN 1845
		15. REGULATARY INFORMATION

## UNITED STATES:

TSCA:

This product is solely for research and development purposes only and may not be used, processed or distributed for a commercial purpose. It may only be handled by technically qualified individuals. Prop 65 Listed Chemicals: No Prop 65 Chemicals. No 313 Chemicals

## CANADA:

DSL/NDSL: Not determined.

COMPONENT DIMETHYL SULFOXIDE WHMIS Classification D2B

### **EUROPEAN UNION:**

PRODUCT RISK PHRASES: None assigned. PRODUCT SAFETY PHRASES: Not applicable. PRODUCT CLASSIFICATION: Not classified

COMPONENT DIMETHYL SULFOXIDE EINECS NUMBER 200-664-3

# 16. OTHER INFORMATION

## Abbreviations

N/A - Data is not applicable or not available SARA - Superfund and Reauthorization Act HMIS - Hazard Material Information System WHMIS - Workplace Hazard Materials Information System NTP - National Toxicology Program OSHA - Occupational Health and Safety Administration IARC - International Agency for Research on Cancer PROP 65 - California Safe Drinking Water and Toxic Enforcement Act of 1986 EINECS - European Inventory of Existing Commercial Chemical Substances

This material is sold for research and development purposes only. It is not for any human or animal therapeutic or clinical diagnostic use. It is not intended for food, drug, household, agricultural, or cosmetic use. An individual technically qualified to handle potentially hazardous chemicals must supervise the use of this material.

The above information was acquired by diligent search and/or investigation and the recommendations are based on prudent application of professional judgment. The information shall not be taken as being all inclusive and is to be used only as a guide. All materials and mixtures may be present unknown hazards and should be used with caution. Since Invitrogen Corporation cannot control the actual methods, volumes, or conditions of use, the Company shall not be held liable for any damages or losses resulting from the handling or from contact with the product as described herein. THE INFORMATION IN THIS MSDS DOES NOT CONSTITUTE A WARRANTY, EXPRESS OR IMPLIED, INCLUDING ANY IMPLIED WARRANTY OF MERCHANTABILITY OR FITNESS FOR ANY PARTCULAR PURPOSE.

# End of Safety Data Sheet