

SECTION 01: PRODUCT AND COMPANY IDENTIFICATION

Product name: Sodium Carbonate
Formula: Na₂CO₃
Chemical family: Carbonates
Synonyms: Disodium carbonate, soda ash and carbonic acid disodium salt.
Product use: For laboratory use only

Manufacturer: CLAISSE
Address: 350, FRANQUET, QUEBEC (QUEBEC) G1P 4P3
CANADA
Phone: +1 418 656-6453
Fax: +1 418 656-1169

Emergency telephone number:
CANUTEC (24h): +1 613 996-6666

SECTION 02: HAZARDS IDENTIFICATION**GHS and (EC) No 1272/2008 classification**

Acute toxicity, Oral (Category 5)
Skin corrosion/irritation (Category 3)
Serious eye damage/Eye irritation (Category 2A).

Label elements:

Pictogram:



Signal word: Warning
SGH 07

Hazard statements:

H303 May be harmful if swallowed.
H316 Causes mild skin irritation.
H319 Causes serious eye irritation.

Precautionary statements:

P261 Avoid breathing dust/fume/gas/mist/vapours/spray.
P264 Wash hands and clothes thoroughly after handling.
P280 Wear protective gloves/protective clothing/eye protection/face protection.
P302+P352 IF ON SKIN: Wash with plenty of soap and water.
P304+P340 IF INHALED: Remove victim to fresh air and keep at rest in a position comfortable for breathing.
P305+P351+P338 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses if present and easy to do. Continue rinsing.

Other hazards

No data available.

SECTION 03: COMPONENT INFORMATION

Compound	Synonym	Molecular formula	Molecular Weight (g/mol)	CAS-No.	EC-No.	Index-No.	Concentration (%)
Sodium Carbonate	See Section 01	Na ₂ CO ₃	105.99	497-19-8	207-838-8	011-005-00-2	100

SECTION 04: FIRST AID MEASURES**Description of first aid measures****General information**

Seek immediate medical advice.
Take affected persons out of danger area and lay down.

After inhalation

In case of unconsciousness place patient stably in side position for transportation.
Supply fresh air. If required, provide artificial respiration. Keep patient warm. If symptoms persist, consult a physician.

After skin contact

Immediately wash with water and soap and rinse thoroughly. If skin irritation persists, consult a physician.

After eye contact

Rinse opened eye for several minutes under running water. If symptoms persist, consult a physician.

After swallowing

Rinse out mouth and then drink plenty of water. If symptoms persist, consult a physician.

Most important symptoms and effects, both acute and delayed

No data available.

Indication of any immediate medical attention and special treatment needed

No data available.

SECTION 05: FIREFIGHTING MEASURES**Suitable extinguishing media**

Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.

Unsuitable extinguishing media

No data available.

Special protective equipment for fire fighters

Wear self-contained breathing apparatus for firefighting if necessary.

Hazardous combustion products

Hazardous decomposition products formed under fire conditions: carbon oxides.

SECTION 06: ACCIDENTAL RELEASE MEASURES**Personal precautions**

Use personal protective equipment. Avoid dust formation. Avoid breathing dust, vapours, fumes or gas. Ensure adequate ventilation. Evacuate personnel to safe areas.

Environmental precautions

Avoid dispersal of spilled material, runoff and contact with soil waterways, drains and sewers.

Methods and materials for containment and cleaning up

Pick up and arrange disposal without creating dust. Sweep up and shovel. Keep in suitable, closed and non-leaking containers for local chemical disposal.

Reference to other sections

See Section 7 for information on safe handling and storage.

See Section 8 for information on exposure controls and personal protection.

See Section 13 for disposal information.

SECTION 07: HANDLING AND STORAGE**Precautions for safe handling**

Provide suction extractors if dust is formed.

Do not inhale dust, smoke or mist.

Avoid contact with the eyes and skin.

Prevent formation of dust.

Prevent formation of aerosols.

Conditions for safe storage, including any incompatibilities

Requirements to be met by storerooms and receptacles: no special measures required.

Information about storage in one common storage facility: not required.

Further information about storage conditions: keep container tightly sealed store receptacle in a well-ventilated area; store in dry conditions.

Specific end use(s)

No data available.

SECTION 08: EXPOSURE CONTROLS/PERSONAL PROTECTION**Control parameters**

No exposure limits have been established.

Exposure Control**Personal protective equipment****General protective and hygienic measures**

Keep away from foodstuffs, beverages and feed. Immediately remove all soiled and contaminated clothing. Wash hands before breaks and at the end of work shifts. Avoid contact with the eyes and skin.

Respiratory protection

For nuisance exposure, use type N95 (US) or type P1 (EU EN 143) particle respirator. For higher level protection, use type OV/AG/P99 (US) or type ABEK-P2 (EU EN 143) respirator cartridges. Use respirators and components tested and approved under appropriate government standards such as NIOSH (US) or CEN (EU).

Hand protection

Handle with gloves. The glove material has to be impermeable and resistant to the product, the substance or preparation. Selection of the glove material must be made considering the penetration times, rates of diffusion and degradation. Gloves must be inspected prior to use. Use proper glove removal technique (without touching glove's outer surface) to avoid skin contact with this product. Dispose of contaminated gloves after use in accordance with applicable laws and good laboratory practices. Wash and dry hands.



Protective gloves.

Eye protection

Wear safety glasses with side shields conforming to EN 166. Use equipment for eye protection tested and approved under appropriate government standards such as NIOSH (US) or EN 166 (EU).



Safety glasses with side shields (EN 166).

Skin and body protection

Wear impervious clothing. The type of protective equipment must be selected according to the concentration and amount of dangerous substance at the specific workplace.

Specific engineering controls

Use mechanical exhaust or laboratory fume hood to avoid exposure.

Environmental exposure controls

Use mechanical exhaust or laboratory fume hood to avoid exposure.

SECTION 09: PHYSICAL AND CHEMICAL PROPERTIES
Appearance

Form: Powder
 Colour: White

Safety data

pH: Basic in solution
 Melting point/freezing point: 851°C/1564°F
 Initial boiling point/boiling range: No data available
 Flash point: No data available
 Flammability: No data available
 Ignition point: No data available
 Autoignition point: No data available
 Lower flammable/explosive limit: No data available
 Upper flammable/explosive limit: No data available
 Vapour pressure: Negligible
 Relative density: 2.53 at 20°C
 Solubility: 215 g/L at 20°C in water
 Partition coefficient n-octanol/water: No data available
 Decomposition temperature: No data available

Viscosity:	No data available
Relative vapour density:	No data available
Odour:	No data available
Odour threshold:	No data available
Evaporation rate:	No data available
Explosive properties:	No data available
Oxidizing properties:	No data available

SECTION 10: STABILITY AND REACTIVITY

Reactivity

The solution in water is alkaline. Decomposes by reaction with strong acids.

Chemical stability

Stable under recommended storage conditions.

Possibility of hazardous reactions

No dangerous reactions known.

Conditions to avoid

Incompatible materials.

Incompatible materials

Finely divided aluminum and strong acids.

Hazardous decomposition products

Hazardous decomposition products formed under fire conditions: carbon oxides and sodium oxides.

SECTION 11: TOXICOLOGICAL INFORMATION

Acute toxicity

Compound	Oral LD50	Inhalation LC50	Dermal LD50	Other
Sodium Carbonate	2800 mg/kg bw	800 mg/m ³ (guinea pig) 1200 mg/m ³ (mice) 2300 mg/m ³ (rat)	>2000 mg/kg bw	No data

Skin corrosion/irritation

Acute skin irritation/corrosion studies were performed with sodium carbonate in accordance with Henkel Hellas S.A. Atalanti. ECETOC, Technical Report No. 66, 1995 (LIT 7835) [IUCLID 2000], EPA 16 CFR 1500.3 and testing comparable to OECD Guideline 404. Based on the results obtained, sodium carbonate has not to be classified and labelled irritating to the skin according to Directive 67/548/EEC (DSD) and Regulation (EC) No 1272/2008 (CLP).

Serious eye damage/eye irritation

Acute eye irritation/corrosion studies were performed with sodium carbonate in accordance with Solvay S.A. Bruxelles Murphy J.C. et al. (1982), Toxicology, 23, p. 281 [IUCLID 2000], EPA 16 CFR 1500.42 and testing comparable to OECD Guideline 405. With respect to eye irritation, sodium carbonate has to be classified into irritating to the eye according to the Regulation (EC) No 1272/2008 (CLP) and according to the Directive 67/548/EEC (DSD).

Respiratory or skin sensitization

No data available.

Germ cell mutagenicity (in vitro) – gene mutation

No data available.

Germ cell mutagenicity (in vivo) – DNA damage and/or repair

No data available.

Carcinogenicity

IARC: No component of this product present at levels greater than or equal to 0.1% is identified as probable, possible or confirmed human carcinogen by IARC.

ACGIH: No component of this product present at levels greater than or equal to 0.1% is identified as a carcinogen or potential carcinogen by ACGIH (A4 group).

Reproductive toxicity

No data available.

STOT – SE (GHS)

No data available.

STOT – RE (GHS)

No data available.

Signs and symptoms of exposure

Burning sensation, cough, wheezing, laryngitis, shortness of breath, headache, nausea and vomiting.

Synergistic effects

No data available.

Additional information

Compound	RTECS
Sodium Carbonate	VZ4050000

SECTION 12: ECOLOGICAL INFORMATION**Toxicity****Aquatic environment toxicity (acute)**

No data available.

Aquatic environment toxicity (long-term)

No data available.

Persistence and degradability

No data available.

Bioaccumulative potential

No data available.

Mobility in soil

No data available.

PBT and vPvB assessment

PBT: Not applicable.

vPvB: Not applicable.

Other adverse effects

Will affect drinking water supplies. The product components are not classified as environmentally hazardous. However, this does not exclude the possibility that large or frequent spills can have harmful or damaging effects on the environment.

SECTION 13: DISPOSAL CONSIDERATIONS**Product disposal**

The generation of waste should be avoided or minimized wherever possible.

Contaminated packaging disposal

Dispose as an unused product.

Waste treatment-relevant information

Disposal of this product, solutions and any by-products should at all times comply with the requirements of environmental protection and waste disposal legislation and any regional and local authority requirements.

Sewage disposal-relevant information

Avoid dispersal of spilled material, runoff and contact with soil, waterways, drains and sewers.

Other disposal recommendations

Contact a licensed professional waste disposal service to dispose of this material.

SECTION 14: TRANSPORT INFORMATION**UN number**

ADR, ADN, IMDG, IATA, TDG, DOT Not applicable.

UN Proper shipping name

Not applicable.

Transport hazard class(es)

ADR, ADN, IMDG, IATA, TDG, DOT Not applicable.

Packing group

ADR/IMDG/IATA/TDG/DOT Not applicable.

Environmental hazards

Environmentally hazardous substance/marine pollutant: No.

Special precaution for user

Not applicable.

Transport in bulk according to Annex II of MARPOL73/78 and the IBC Code

Not applicable.

SECTION 15: REGULATORY INFORMATION**Safety, health and environmental regulations/legislation specific for the substance or mixture**

This safety data sheet complies with the requirements of regulation (EC) No. 1907/2006.

Chemical safety assessment

A chemical safety assessment has not been carried out.

DSL status

All components of this product are on the Canadian DSL list.

WHMIS classification

D2B Toxic material Moderate eye irritant

This product has been classified in accordance with the hazard criteria of the CPR and the MSDS contains all the information required by the CPR.

HMIS classification

Health hazard:	1
Chronic health hazard:	*
Flammability:	0
Physical hazards:	0

Potential health effects

Inhalation:	May cause respiratory irritation.
Skin:	May cause skin irritation.
Eyes:	Causes serious eye irritation.
Ingestion:	May be harmful if swallowed.

SECTION 16: OTHER INFORMATION**Date of issue**

2016-08-15

Notice to the reader

To the best of our knowledge, the information contained herein is accurate. However, neither the above named supplier nor any of its subsidiaries assumes any liability whatsoever for the accuracy or completeness of the information contained herein. Final determination of suitability of any material is the only responsibility of the user. All materials may present unknown hazards and should be used with caution. Although certain hazards are described herein, we cannot guarantee that these are the only.

Abbreviations and acronyms

ACGIH:	American Conference of Governmental Industrial Hygienists
ADR:	European Agreement concerning the International Carriage of Dangerous Goods by Road
ADN:	European Agreement concerning the International Carriage of Dangerous Goods by Inland Waterway
b.w.:	Body weight
CAA:	Clean Air Act
CAS:	Chemical Abstracts Service (division of the American Chemical Society)
CEN:	European Committee for Standardization
CERCLA:	Comprehensive Environmental Response, Compensation and Liability Act
CFR:	Code of Federal Regulations
CLP:	Classification, Labelling and Packaging
CPR:	Controlled Products Regulations
DNEL:	Derived No-Effect Level
DOT:	Department of Transportation
DSL:	Domestic Substance List
EINECS:	European Inventory of Existing Commercial Chemical Substances
GHS:	Globally Harmonized System of Classification and Labelling of Chemicals
HDPE:	High Density PolyEthylene
HEPA:	High Efficiency Particulate Air
HMIS:	Hazardous Material Information System
IARC:	International Agency for Research on Cancer
IATA:	International Air Transport Association

IBC:	Intermediate Bulk Container
IDLH:	Immediately Dangerous to Life or Health Concentrations
IMDG:	International Maritime Dangerous Goods Code
LC50:	Median Lethal Concentration
LD50:	Median Lethal Dose
LOAEL:	Lowest Observed Adverse Effect Level
LOEC:	Lowest Observable Effect Concentration
MARPOL:	MARine POLLution
NIOSH:	The National Institute for Occupational Safety and Health
NOAEL:	No Observed Adverse Effect Level
NOEC:	No Observable Effect Concentration
OECD:	Organisation for Economic Co-operation and Development
OSHA:	Occupational Safety and Health Administration
PBT:	Persistent Bioaccumulative and Toxic
PEL:	Permissible Exposure Limits
PNEC:	Predicted No-Effect Concentration
RCF:	Refractory Ceramic Fibers
RTECS:	Registry of Toxic Effects of Chemical Substances
SARA:	Superfund Amendments and Reauthorization Act
SCP:	Standards Completion Program (NIOSH/OSHA)
STEL:	Short Term Exposure Limit
STOT – RE:	Specific Target Organ Toxicity – Repeated exposure
STOT – SE:	Specific Target Organ Toxicity – Single exposure
TDG:	Transport of Dangerous Goods
TLV:	Threshold Limit Value
TSCA:	Toxic Substances Control Act
TWA:	Time-Weighted Average exposure value
UN:	United Nations
vPvB:	very Persistent and very Bioaccumulative
WHMIS:	Workplace Hazardous Materials Information System