

SECTION 01: PRODUCT AND COMPANY IDENTIFICATION

Product name: Sodium Iodide
Formula: NaI
Chemical family: Iodide
Synonyms: No data available
Recommended use: Non wetting agent for sample preparation by fusion
Restrictions on 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, WHMIS 2015 and (EC) No 1272/2008 classification

Skin irritation (Category 2)
 Eye irritation (Category 2A)
 Acute aquatic toxicity (category 1)
 STOT RE (Category 1)

Label elements:

Pictograms:



Signal word:

Health hazard
(GHS08)

Environmental hazard
(GHS09)

Hazard statements:

H315	Causes skin irritation.
H319	Causes serious eye irritation.
H372	Causes damage to thyroid gland through prolonged or repeated exposure.
H400	Very toxic to aquatic life.

Precautionary statements:

P260	Do not breathe dust/fume/gas/mist/vapours/spray.
P264	Wash hands thoroughly after handling.
P270	Do not eat, drink or smoke when using this product.
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

Compounds	Molecular formula	Synonym	Molecular Weight (g/mol)	CAS-No.	EC-No.	Concentration (%)
Sodium iodide	Nal	-	149.89	7681-82-5	231-679-3	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.

Innapropriate extinguishing media

No data available.

Special protective equipment and precaution for fire fighters

Wear self-contained breathing apparatus for firefighting if necessary.

Special hazards arising from the substance or mixture

Hazardous decomposition products formed under fire conditions: sodium oxide and hydrogen iodide.

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.

See Section 8 for information on personal protective equipment.

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

Compound	Exposure Limits	Basis	Entity
Sodium iodide	0.01 mg/m ³	TWA	ACGIH

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 P95 (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 engineering controls

No data available.

SECTION 09: PHYSICAL AND CHEMICAL PROPERTIES
Appearance

Form: Powder
 Colour: White

Safety data

Odour: No data available
 Odour threshold: No data available
 pH: No data available
 Melting point/freezing point: 661°C/1222°F
 Initial boiling point/boiling range: 1304°C/2379°F
 Flash point: No data available
 Evaporation rate: No data available
 Flammability: No data available
 Lower flammable/explosive limit: No data available
 Upper flammable/explosive limit: No data available
 Vapour pressure: No data available
 Relative density: No data available
 Solubility: No data available
 Partition coefficient n-octanol/water: No data available

Autoignition point:	No data available
Ignition point:	No data available
Decomposition temperature:	No data available
Viscosity:	No data available
Relative vapour density:	No data available
Explosive properties:	No data available
Oxidizing properties:	No data available

SECTION 10: STABILITY AND REACTIVITY

Reactivity

No data available.

Chemical stability

Stable under recommended storage conditions.

Possibility of hazardous reactions

No dangerous reactions known.

Conditions to avoid

Exposure to light and air may affect the quality of the product.

Incompatible materials

No data available.

Hazardous decomposition products

Hazardous decomposition products formed under fire conditions: sodium oxide and hydrogen iodide.

SECTION 11: TOXICOLOGICAL INFORMATION

Acute toxicity

Compounds	Oral LD ₅₀	Inhalation LC ₅₀	Dermal LD ₅₀	Other
Sodium Iodide	4340 mg/kg, rat	No data	No data	No data

Repeated exposure toxicity

Compounds	Oral DNEL	Inhalation DNEL	Dermal DNEL	Other
Sodium Iodide	0.01 mg/kg bw/day	No data	No data	No data

Skin corrosion/irritation

No data available.

Serious eye damage/eye irritation

No data available.

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.

Reproductive toxicity

No data available.

STOT – SE (GHS)

No data available.

STOT – RE (GHS)

No data available.

Potential health effects

Inhalation: May be harmful if inhaled. Causes respiratory tract irritation.

Ingestion: May be harmful if swallowed.

Skin: May be harmful if absorbed through skin. Causes skin irritation.

Eyes: Causes eye irritation.

Signs and symptoms of exposure

Prolonged exposure to iodides may produce iodism in sensitive individuals. Symptoms of exposure include: skin rash, running nose, headache and irritation of the mucous membrane. For severe cases the skin may show pimples, boils, hives, blisters and black and blue spots. Iodides are readily diffused across the placenta. Neonatal deaths from respiratory distress secondary to goiter have been reported. Iodides have been known to cause drug-induced fevers, which are usually of short duration. To the best of our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated.

Synergistic effects

No data available.

Additional information

Compounds	RTECS
Sodium Iodide	WB6475000

SECTION 12: ECOLOGICAL INFORMATION

Toxicity

Aquatic toxicity

Compounds	NOEC	LOEC
Sodium Iodide	No data	No data

Persistence and degradability

No data available.

Bioaccumulative potential

No data available.

Mobility in soil

Water hazard class 1 (German regulation, self-assessment): slightly hazardous for water. Do not allow undiluted product or large quantities of it to reach ground water, water course or sewage system.

PBT and vPvB assessment

PBT: Not applicable.
vPvB: Not applicable.

Other adverse effects

Will affect drinking water supplies. The product components is classified as environmentally hazardous.

SECTION 13: DISPOSAL CONSIDERATIONS**Product disposal**

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

Do not disperse or spill the material. Runoff and contact with soil, waterways, drains and sewers must not happen. Disposal of this solutions should at all times comply with the requirements of environmental protection and waste disposal legislation and any regional and local authority requirements.

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, DOT, TDG UN3077

UN Proper shipping name

ENVIRONMENTALLY HAZARDOUS SUBSTANCE, SOLID, N.O.S. (Sodium iodide)

Transport hazard class(es)

ADR, ADN, IMDG, IATA, DOT, TDG 9

Packing group

ADR, ADN, IMDG, IATA, DOT, TDG III

Environmental hazards

Environmentally hazardous substance/marine pollutant: Yes

Special precaution for user

EHS-Mark required (ADR 2.2.9.1.10, IMDG code 2.10.3) for single packagings and combination packagings containing inner packagings with Dangerous Goods > 5L for liquids or > 5kg for solids.

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

Not applicable.

SECTION 15: REGULATORY INFORMATION**Emergency overview**

Target organs: Thyroid gland, blood, bone marrow.

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 1988

D2B Toxic material causing other toxic effect Moderate skin irritant
Moderate eye irritant

HMIS 1988 classification

Health hazard: 2
Chronic health hazard: *
Flammability: 0
Physical hazards: 0

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

2017-02-14

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
OEL:	Occupational Exposure Level
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