

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.
cleaning@alconox.com

- alconox.com
- distributors
 - technical information
 - samples
 - new developments
 - white papers
 - prices

LUMINOX[®]

Low-Foaming Neutral Cleaner

- Low-foaming for use in mechanical and spray washers
- Neutral pH for non-corrosive cleaning, safe handling and easy waste disposal
- Concentrated for economical use
- Free rinsing to eliminate interfering residues
- Biodegradable and phosphate free

Used to clean: Aluminum and soft metal surfaces, medical instruments in aluminum trays, alkaline-sensitive lab and glassware, sensitive aerospace materials, pharmaceutical process equipment, medical devices, electronic components.

Used to remove: Oils, dirt, grime, adhesives and uncured epoxy, particulates, residues, fluxes, chemical solvents, machining oils.

Surfaces cleaned: Corrosion inhibited formulation recommended for glass, ferrous metals, non-ferrous metals, aluminum, copper, zinc, and other soft metals, ceramic, plastic, circuit boards and electronic components.

Cleaning method: Soak tanks, manual cleaning, ultrasonic tanks, mechanical washers, washer/sanitizers, spray washers, CIP systems.

Directions: For manual, soak and ultrasonic use, make a 1 – 2% solution (3 – 5 Tbsp/gal, 1 1/4 – 2 1/2 oz/gal or 10 – 20 ml/L) in warm or hot water. Protective gloves and eyewear are recommended. For spray machines, add detergent at the hot water wash cycle to make a 2% solution (2 1/2 oz per gal, or 20 ml per L). For difficult soils raise water temperature and use more detergent. RINSE THOROUGHLY—preferably with running water. For critical cleaning do final or all rinsing

in distilled, deionized or purified water. Corrosion testing may be advisable. Material compatibility guidelines available on request.

Convenient Sizes	Alconox Cat. #
Case of 4 x 1 gal (4 x 3.8L)	1901
5 gal (19L) jerrycan	1905
15 gal (57L) drum	1915
55 gal (208L) drum	1955
1 gal makes 20–50 gal (1 L makes 20–50L) solution	



LUMINOX passes inhibitory residue test for water analysis and is FDA certified. LUMINOX is available from leading laboratory, hospital, clinical and industrial suppliers. To find a distributor for Alconox Inc. detergents, use the Locate Dealers search tool at alconox.com. To request samples, use the Order Samples form at alconox.com. Or call Alconox Inc. at +914.948.4040, today.

LUTB.4

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.

cleaning@alconox.com

alconox.com

- distributors
- technical information
- samples
- new developments
- white papers
- prices

Physical Data	Typical Value
pH (as is)	7.0
Surface Tension 1% Sol'n (dyne/cm)	34
Conductivity 1% Sol'n (mS/cm at 21 degrees C)	0.4
Form	Liquid
Color	Clear
Solubility in Water	Completely soluble in all proportions
Hard Water Effectiveness	Highly effective
Biodegradability	Yes, meets EC648/2004
Foam Tendency	Low foaming
Shelf Life	42 months from end of month of manufacture
Fragrance and Dye Content	0%
Phosphate Content (as phosphorus)	0%
Flash Point (degrees F)	>200 degrees F
Density (lb/gal)	8.75

Chemical Description

Luminox is an aqueous, low-ionic content, phosphate-free blend of organic solvents, mixture of nonionic and anionic surfactants, and rinse aids.

Cleaning Validation and Residue Detection

To quantify trace levels of Luminox for cleaning validation purposes, use TOC, or HPLC. For details see the Cleaning Validation page in Support at alconox.com. For lab accreditation tests to show the detergent is rinsed away, see the Accreditation Residue Detection in the Support section of alconox.com.

Health and Safety Information

OSHA Hazardous Ingredients	None
RCRA Hazard Class	Nonhazardous
EPA Priority Pollutants	None
DOT Hazard Class	Nonhazardous
Flammability	Nonflammable
Latex Content	None in product or packaging.
Irritation	Cat. 2 skin, Cat. 2A eye irritant
Oral Toxicity	LD ₅₀ > 5000 mg/kg oral rat.
Inhalation Toxicity	Nonirritating
VOC Content	Less than 50 g/liter even at 10% use concentration.
GMO Content	None
BSE/TSE Source	None
2007/68/EC Annex IIIa Allergens	None
REACH Art. 57 (SVHC)	None
RoHS	No RoHS hazardous ingredients.
Carcinogenicity	NTP = No IARC = No OSHA = No
Ingredients on TSCA Inventory	Yes
EC 907/2006	Contains no enzymes, optical brighteners, perfumes, allergenic fragrances, or preservative agents.

Precautions

Avoid contact with skin, eyes and clothing. Do not take internally. In case of contact immediately flush skin with plenty of water. For eyes, flush with plenty of water for at least 15 minutes and get medical attention immediately. A Safety Data Sheet is available at alconox.com. Contact Alconox Inc. for purchase specifications. Information presented is typical. Not to be taken as specifications. Typical data is not a specification.

While the information in this report should not be considered to be a product warranty, we urge you to investigate, test and verify the suitability of Alconox Inc. detergents for your specific application. We, of course, can not give permission to use, or recommend the use of, our detergents where they infringe patents. No representation or warranty is made as to the safety of products or materials mentioned under the Federal Food Additives Amendment of 1958.

LUTB.4