

Globally Harmonized System Pictogram Reference Table

The following summarizes new label elements per the GHS classification guidelines. Each summary describes the previous label element compared to the new, and may include examples of the new elements.

Pictogram

The pictogram, previously known as a “symbol”, has a black symbol on a white background with red border and is in shape of a square set at a point or a diamond. Some of the lower hazard categories may not have an assigned pictogram. For health hazard precedence may apply for example “Exclamation mark” pictogram may not appear for skin or eye irritation if the “Corrosive” pictogram applies. GHS has a total of nine pictograms.

Signal words

Signal word is a new implementation of GHS and indicates the level of hazard severity. The signal words “Danger” or “Warning” are assigned to a specific hazard class. In case of multiple hazard classes for one product, the word “Danger” takes precedence over “Warning”. Some of the lower hazard categories may not be assigned a signal word.

Hazard statements

The hazard (**H**) statement, previously known as a “Risk Phrase”, is assigned to each hazard class and category. The GHS consists of 3 hazard groups: physical, health, and environmental. Each group has multiple hazard classes, further containing categories and sub categories. The H statement has a unique code consisting of the letter “H”, a number for the hazard group followed by two sequential numbers. The assigned hazard group numbers are:

- 2 – Physical hazard
- 3 – Health hazard
- 4 – Environmental hazard





The examples of **H** statement codes are:

- 2 – Physical hazard - **H226** “Flammable liquid and vapor”
- 3 – Health hazard - **H303** “May be harmful if swallowed”
- 4 – Environmental hazard - **H401** “Toxic to aquatic life”







For your reference, the table within the following pages outlines the various combinations of Hazard Class, Pictograms, Signal Words, Hazard Statement Codes, and Description of Hazard Statements in relation to each hazard group. Further information can be found by following the reference web link at the bottom of each page.

Globally Harmonized System Pictogram Reference Table






GHS Health Hazards

Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Acute Toxicity Oral Category 1		Danger	H300	Fatal if swallowed
Acute Toxicity Dermal Category 1			H310	Fatal in contact with skin
Acute Toxicity Inhalation Category 1			H330	Fatal if inhaled
Acute Toxicity Oral Category 2			H300	Fatal if swallowed
Acute Toxicity Dermal Category 2			H310	Fatal in contact with skin
Acute Toxicity Inhalation Category 2			H330	Fatal if inhaled
Acute Toxicity Oral Category 3			H301	Toxic if swallowed
Acute Toxicity Dermal Category 3			H311	Toxic in contact with skin
Acute Toxicity Inhalation Category 3			H331	Toxic if inhaled
Acute Toxicity Oral Category 4		Warning	H302	Harmful if swallowed
Acute Toxicity Dermal Category 4			H312	Harmful in contact with skin
Acute Toxicity Inhalation Category 4			H332	Harmful if inhaled
Acute Toxicity Oral Category 5	None	Warning	H303	May be harmful if swallowed
Acute Toxicity Dermal Category 5			H313	May be harmful In contact with skin
Acute Toxicity Inhalation Category 5			H333	May be harmful if inhaled
Skin Corrosion /Irritation Category 1		Danger	H314	Causes severe skin burns and eye damage
Skin Corrosion /Irritation Category 2		Warning	H315	Causes skin irritation
Skin Corrosion /Irritation Category 3	None	Warning	H316	Causes mild skin irritation

Globally Harmonized System Pictogram Reference Table




Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Serious eye damage/eye irritation Category 1		Danger	H318	Causes serious eye damage
Serious eye damage/eye irritation Category 2/2A		Warning	H319	Causes serious eye irritation
Serious eye damage/eye irritation Category 2B	None	Warning	H320	Causes eye irritation
Respiratory Sensitization Category 1		Danger	H334	May cause allergy or asthma symptoms or breathing difficulties if inhaled
Respiratory Sensitization Category 1A				
Respiratory Sensitization Category 1B				
Skin Sensitization Category 1		Warning	H317	May cause an allergic skin reaction
Skin Sensitization Category 1A				
Skin Sensitization Category 1B				
Germ cell mutagenicity Category 1		Danger	H340	May cause genetic defects
Germ cell mutagenicity Category 2		Warning	H341	Suspected of causing genetic defects
Carcinogenicity Category 1		Danger	H350	May cause cancer
Carcinogenicity Category 2		Warning	H351	Suspected of causing cancer

Globally Harmonized System Pictogram Reference Table

Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Reproductive Toxicity Category 1		Danger	H360	May damage fertility or the unborn child
Reproductive Toxicity Category 2		Warning	H361	Suspected of damaging fertility or the unborn child
Reproductive Toxicity-Effect via lactation	None	None	H362	May cause harm to breast-fed children
Specific Target Organ toxicity - single exposure Category 1		Danger	H370	Causes damage to organs
Specific Target Organ toxicity - single exposure Category 2		Warning	H371	May cause damage to organs
Specific Target Organ toxicity - single exposure Category 3		Warning	H335 or H336	May cause respiratory irritation Or May cause drowsiness or dizziness
Specific Target Organ toxicity - repeated exposure Category 1		Danger	H372	Causes damage to organs through prolonged or repeated exposure
Specific Target Organ toxicity - repeated exposure Category 2		Warning	H373	May cause damage to organs through prolonged or repeated exposure
Aspiration Hazard Category 1		Danger	H304	May be fatal if swallowed and enters airways
Aspiration Hazard Category 2		Warning	H305	May be harmful if swallowed and enters airways

Globally Harmonized System Pictogram Reference Table

GHS Environmental Hazards









Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Hazardous to aquatic environment Short term /Acute Category 1		Warning	H400	Very toxic to aquatic life
Hazardous to aquatic environment Short term /Acute Category 2	None	None	H401	Toxic to aquatic life
Hazardous to aquatic environment Short term /Acute Category 3	None	None	H402	Harmful to aquatic life
Hazardous to aquatic environment Long term/Chronic Category 1		Warning	H410	Very toxic to aquatic life with long lasting effects
Hazardous to aquatic environment Long term/Chronic Category 2		None	H411	Toxic to aquatic life with long lasting effects
Hazardous to aquatic environment Long term/Chronic Category 3	None	None	H412	Harmful to aquatic life with long lasting effects
Hazardous to aquatic environment Long term/Chronic Category 4	None	None	H413	May cause long lasting harmful effects to aquatic life
Hazardous to ozone layer Category 1		Warning	H420	Harms public health and the environment by destroying ozone in the upper atmosphere

Globally Harmonized System Pictogram Reference Table








GHS Physical Hazard

Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Unstable explosive		Danger	H200	Unstable explosives
Explosives Division 1.1			H201	Explosive; mass explosion hazard
Explosives Division 1.2			H202	Explosive, severe projection hazard
Explosives Division 1.3			H203	Explosive; fire, blast or projection hazard
Explosives Division 1.4		Warning	H204	Fire or projection hazard
Explosives Division 1.5	None	Danger	H205	May mass explode in fire
Explosives Division 1.6	None	None	None	None
Flammable Gases Category 1		Danger	H220	Extremely flammable gas
Flammable Gases Category 2	None	Warning	H221	Flammable gas
Aerosols Category 1		Danger	H222, H229	Extremely flammable aerosol; Pressurized container: may burst if heated
Aerosols Category 2		Warning	H223, H229	Flammable aerosol; Pressurized container: may burst if heated
Aerosols Category 3	None	Warning	H229	Pressurized container: may burst if heated
Oxidizing gases Category 1		Danger	H270	May cause or intensify fire; oxidiser
Gases under pressure - Compressed gas		Warning	H280	Contains gas under pressure; may explode if heated
Gases under pressure - Liquefied gas				
Gases under pressure - Dissolved gas				
Gases under pressure - Refrigerated gas			H281	Contains refrigerated gas; may cause cryogenic burns or injury
Flammable Liquids Category 1		Danger	H224	Extremely flammable liquid and vapour
Flammable Liquids Category 2		Danger	H225	Highly flammable liquid and vapour
Flammable Liquids Category 3		Warning	H226	Flammable liquid and vapour
Flammable Liquids Category 4	None	Warning	H227	Combustible Liquid

Globally Harmonized System Pictogram Reference Table

Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Flammable Solids Category 1		Danger	H228	Flammable solid
Flammable Solids Category 2		Warning		
Self-reactive substances and mixtures Type A		Danger	H240	Heating may cause an explosion
Self-reactive substances and mixtures Type B	 	Danger	H241	Heating may cause a fire or explosion
Self-reactive substances and mixtures Type C and D		Danger	H242	Heating may cause a fire
Self-reactive substances and mixtures Type E and F		Warning		
Self-reactive substances and mixtures Type G	None	None	None	None
Pyrophoric liquids Category 1		Danger	H250	Catch fire spontaneously if exposed to air
Pyrophoric solids Category 1				
Self-heating substances and mixtures Category 1		Danger	H251	Self-heating; may catch fire
Self-heating substances and mixtures Category 2		Warning	H252	Self-heating in large quantities; may catch fire
Substances and mixtures which in contact with water emit flammable gases Category 1		Danger	H260	In contact with water releases flammable gases which may ignite spontaneously
Substances and mixtures which in contact with water emit flammable gases Category 2			H261	In contact with water releases flammable gases
Substances and mixtures which in contact with water emit flammable gases Category 3		Warning	H261	In contact with water releases flammable gases

Globally Harmonized System Pictogram Reference Table

Hazard Class and Categories	Pictogram	Signal Word	Hazard Statement Code	Description of Hazard Statement
Oxidizing liquids Category 1		Danger	H271	May cause fire or explosion; strong oxidizer
Oxidizing liquids Category 2		Danger	H272	May intensify fire; oxidizer
Oxidizing liquids Category 3		Warning		
Oxidizing Solids Category 1		Danger	H271	May cause fire or explosion; strong oxidizer
Oxidizing Solids Category 2		Danger	H272	May intensify fire; oxidizer
Oxidizing Solids Category 3		Warning		
Organic peroxides Type A		Danger	H240	Heating may cause an explosion
Organic peroxides Type B	 	Danger	H241	Heating may cause a fire or an explosion
Organic peroxides Type C and D		Danger	H242	Heating may cause fire
Organic peroxides Type E and F		Warning	H242	Heating may cause fire
Organic peroxides Type G	None	None	None	No hazard statement
Corrosive to metal Category 1		Warning	H290	May be corrosive to metal