



SOLUBILITY RULES CHART

	GROUP I ALKALI METALS				GROUP II ALKALINE EARTH METALS			TRANSITION METALS					POST-TRANSITION METALS	
	Ammonium (NH ₄ ⁺)	Lithium (Li ⁺)	Sodium (Na ⁺)	Potassium (K ⁺)	Magnesium (Mg ²⁺)	Calcium (Ca ²⁺)	Barium (Ba ²⁺)	Iron (II) (Fe ²⁺)	Iron (III) (Fe ³⁺)	Copper (II) (Cu ²⁺)	Silver (Ag ⁺)	Zinc (Zn ²⁺)	Lead (II) (Pb ²⁺)	Aluminum (Al ³⁺)
Fluoride (F ⁻)	SOLUBLE	SLIGHTLY SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	INSOLUBLE	SLIGHTLY SOLUBLE	SLIGHTLY SOLUBLE	SLIGHTLY SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	SLIGHTLY SOLUBLE
Chloride (Cl ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	SOLUBLE	INSOLUBLE	SOLUBLE
Bromide (Br ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	SOLUBLE	SLIGHTLY SOLUBLE	SOLUBLE
Iodide (I ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE			INSOLUBLE		INSOLUBLE	SOLUBLE
Chlorate (ClO ₃ ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE		SOLUBLE		SOLUBLE		SOLUBLE	SOLUBLE
Hydroxide (OH ⁻)		SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	SLIGHTLY SOLUBLE	SOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	SLIGHTLY SOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE
Sulfite (SO ₃ ²⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	INSOLUBLE				INSOLUBLE	INSOLUBLE	INSOLUBLE	
Sulfate (SO ₄ ²⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SLIGHTLY SOLUBLE	INSOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SLIGHTLY SOLUBLE	SOLUBLE	INSOLUBLE	SOLUBLE
Carbonate (CO ₃ ²⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE		INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	
Nitrite (NO ₂ ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE				INSOLUBLE		SOLUBLE	
Nitrate (NO ₃ ⁻)	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE	SOLUBLE
Phosphate (PO ₄ ³⁻)	SOLUBLE	INSOLUBLE	SOLUBLE	SOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE	INSOLUBLE