

9 Fast Facts About Silver

Atomic Number
47

From the Anglo-Saxon word "siolfur." The symbol Ag comes from the Latin "argentum." – Both mean "silver."

Has been known since ancient times

Can be used in:

- Jewelry and Coins
- Electronics
- Photography
- Medical Industry
- Plastics

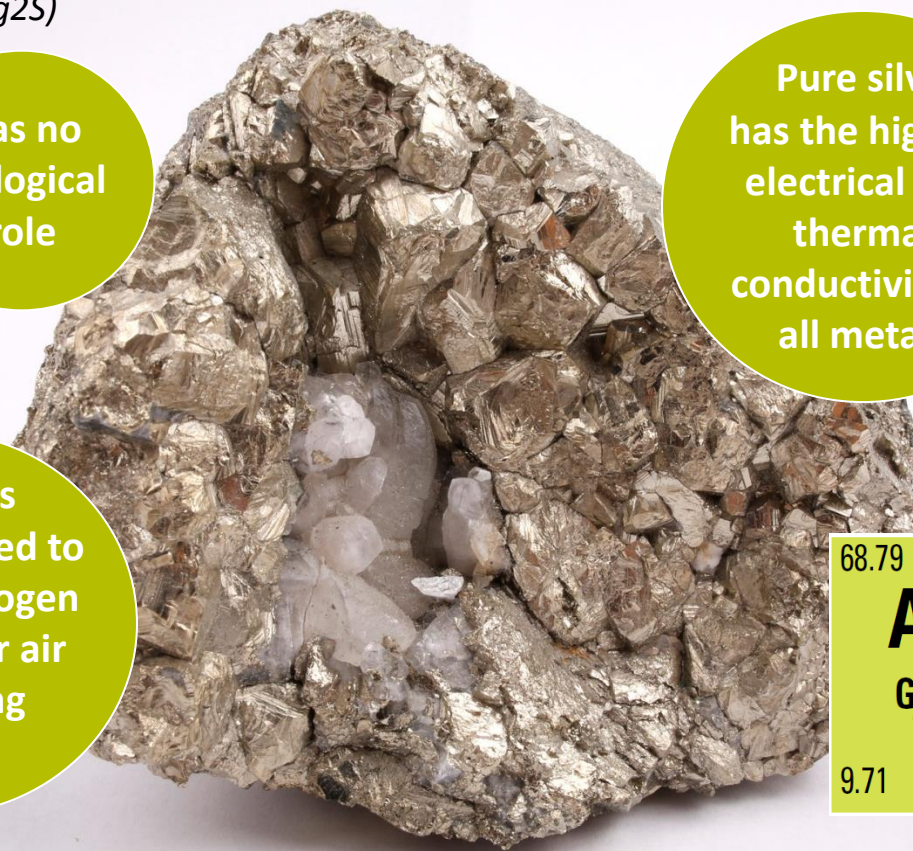
Has an alchemical symbol

Silver is found sometimes as a free metal as well as in ores such as argentite (silver sulphide, Ag₂S)

Has no biological role

Pure silver has the highest electrical and thermal conductivity of all metals

Tarnishes when exposed to ozone, hydrogen sulphide, or air containing sulphur



68.79	77.97
Au	
Gold	
79	
9.71	11.44

Sterling silver is the standard of quality for articles containing 92.5% silver and 7.5% copper (and/or other alloys).

References:
www.webelements.com/silver
www.usgs.gov

©2017 Thermo Fisher Scientific Inc.
All rights reserved.

Delivering Value to the Mining Industry

www.thermofisher.com/mining

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific