

PRODUCT INFORMATION



Heavy Metals Mixture 5 (CRM)

Item No. 41850

Supplied as: A solution in 5% nitric acid (100 µg/ml each compound)

Fill volume: 30 ml

Storage: Room temperature

Stability: ≥3 years

Special conditions: Vortex or sonicate before use

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Heavy Metals Mixture 5 (CRM) (Item No. 41850) is a certified reference material mixture containing heavy metals, which have been found as environmental contaminants in water, soil, plants, and *Cannabis* products.¹⁻⁷ The mixture contains arsenic (As), cadmium (Cd), lead (Pb), mercury (Hg), and chromium (Cr). Levels of these heavy metals are frequently monitored in *Cannabis* samples and have been monitored in *C. sativa* plants used in phytoremediation of cadmium-polluted soils.⁸⁻¹² The mixture is supplied in a plastic bottle with a 30 ml fill volume. It is designed for use in atomic absorption, inductively coupled plasma (ICP), and ICP-MS applications for the detection and quantification of these heavy metals in environmental monitoring and action limit testing. This product is intended for research and environmental testing applications.

This mixture contains five heavy metals found as environmental contaminants in water, soil, plants, and *Cannabis* products. Included are:

Item No. 41850	Heavy Metals Mixture 5 (CRM)
Item Name	Concentration (µg/ml)
Arsenic	100
Cadmium	100
Lead	100
Mercury	100
Chromium	100

References

1. Durum, W.H. and Hem, J.D. *Ann. N.Y. Acad. Sci.* **199**, 26-36 (1972).
2. Williams, C. *J. Sci. Food Agric.* **27(6)**, 561-570 (1976).
3. Laza, A., Orozco, E., and Fernández-Baldo, M.A. *Microchem. J.* **158**, 105265 (2020).
4. Varun, M., D'Souza, R., Pratas, J., et al. *Environ. Sci. Pollut. Res. Int.* **19(1)**, 269-281 (2012).
5. Moir, D., Rickert, W.S., Levasseur, G., et al. *Chem. Res. Toxicol.* **21(2)**, 494-502 (2008).
6. Gardener, H., Wallin, C., and Bowen, J.A. *Sci. Total Environ.* **851(Pt. 1)**, 158110 (2022).
7. Boszke, L., Kowalski, A., Astel, A., et al. *Earth Sci.* **55(5)**, 1075-1087 (2008).
8. Craven, C.B., Wawryk, N., Jiang, P., et al. *J. Environ. Sci. (China)* **85**, 82-93 (2019).
9. Sharma, P., Tripathi, S., Purchase, D., et al. *J. Environ. Chem. Eng.* **9(4)**, 105547 (2021).
10. Kos, B. and Leštan, D. *Plant and Soil* **263**, 43-51 (2004).
11. Goldman, S., Bramante, J., Vrdoljak, G., et al. *J. Liq. Chromatogr.* **44(9-10)**, 403-420 (2021).
12. Linger, P., Müssig, J., Fischer, H., et al. *Ind. Crops Prod.* **16(1)**, 33-42 (2002).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/28/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM