



Product Description: ZeroC™ Dried Plant Matrix

ZeroC is a novel cannabinoid-free *Cannabis* reference matrix designed to meet the rigorous demands of laboratories, researchers, and government agencies. It serves as an essential tool in the field of *Cannabis* science for quality control, method validation, and accurate analyte quantification.

ZeroC allows laboratories to reliably check the effectiveness of their analytical instruments, validate their test methods, and combat the problem of "laboratory shopping" by establishing a standardized blank matrix. This *Cannabis* plant-based reference material serves as a comprehensive solution for investigating common interferences associated with testing cannabinoids, pesticides, mycotoxins, heavy metals, terpenes, residual solvents, and microbials.

ZeroC is a naturally produced and non-extracted form of *Cannabis*, void of any major or minor phytocannabinoids including THC, CBN, CBG, CBD, and THCV. Additionally, it is deficient in many common *Cannabis* monoterpenes and sesquiterpenes. Each lot of reference matrix material is thoroughly characterized and includes certificates of analysis, offering insights into the composition, purity, and traceability.

ZeroC empowers labs to adhere to regulatory compliance and establish robust quality assurance programs that exceed industry benchmarks.

WARNING: ZeroC is intended for laboratory use only and should be handled by trained professionals following appropriate safety protocols. This product is not for human or veterinary use.