

Title: **Workflow solution for steroids and antidoping analysis in urine with GC-MS/MS and GC-HRAM.**

Authors: Petra Gerhards 1, Dr. Daniela Cavagnino 2, Inge de Dobbeleer 3.

1 Thermo Fisher Scientific, Dreieich, Germany, 2 Thermo Fisher Scientific, Milano, Italy, 3 Thermo Fisher Scientific, Breda, The Netherlands

Presenter: Daniela Cavagnino

Purpose

The detection of performance enhancing drugs in sport competition is a growing analytical challenge due to the ongoing development of new drugs and the continuous abuse of older and discontinued drugs found in the black market.

This presentation will show a complete workflow solution including sample preparation for different doping substances as well as confirmation and quantification using GC-MS/MS and GC- HRAM.

Methods and materials

For sample preparation, a C18 and a cation exchange mechanism on a SPE (Solid Phase Extraction) cartridge is described explaining the advantages of SPE compared to LLE (Liquid Extraction).

For analysis, results from a GC coupled with a triple quadrupole MS are shown and to perform screening and confirmation in one analysis, data are shown on a GC-HRAM.

A summary of results

SPE yields cleaner extracts and can offer a concentration factor of up to 120 times, gaining better detection limits.

GC-MS/MS offers good selectivity in the urine and results on detection limits and robustness are shown obtained by the Rio de Janeiro Doping control laboratory.

For the non-targeted approach for new substances results are shown on a GC-HRAM within 1ppm mass accuracy across a broad linearity range, including good sensitivity, and linearity for a large number of compounds.

Statement of conclusions reached

The method shows detection limits below the minimum required performance limits for both a targeted approach and non-targeted approach with various GCMS technologies. Especially for new psychoactive substances.