



Diagnosing EASY-Spray chromatographic problems

Authors

CCS Center of Excellence (COE)
Application Scientists
Thermo Fisher Scientific

Keywords

EASY-Spray, nanoLC, proteomics,
peptides

Introduction

Chromatographic problems with Thermo Scientific™ EASY-Spray™ columns are best evaluated using a simple diagnostic standard such as Thermo Scientific™ Pierce™ Peptide Retention Time Calibration mixture (PRTC), Thermo Scientific™ Pierce™ BSA (bovine serum albumin) Protein Digest Standard, or Thermo Scientific™ Dionex™ Cytochrome C digest.

This document describes the steps to collect and analyze the appropriate data for diagnosing issues with EASY-Spray columns to facilitate in the identification of system contamination, sample quality issues and overall column performance.

Use of diagnostic standards for EASY-Spray columns

1. Diagnostic standards

- Use a diagnostic standard that allows the extraction of a known peptide m/z
- The standard used in EASY-Spray quality testing is Thermo Scientific™ Pierce™ Bovine Serum Albumin – Tryptic Digest, MS grade (PN 88341)
- Alternate standards that may be used include Pierce PRTC mixture (PN 88320) and Cytochrome C digest (PN 161089)

2. Instrument Settings

- If using the BSA standard, use the LC-MS Analysis Parameters on the Quality Assurance Report shipped with the column as your reference
- For example, column ES804A LC settings:

Mobile phase A:	100% Water, 0.1% Formic acid	Mobile phase B:	100% Acetonitrile, 0.1% Formic acid
Flow rate:	300 nL/min	Column temperature:	35 °C
Gradient:	5–40% B in 15 min 40–95% B in 5 min 95% B for 5 min	Run time:	25 min
Equilibration:	5% B for 20 min		

- MS settings

Acquisition range:	400–2000 <i>m/z</i>	Spray voltage:	1.9 kV
Scan rate:	Normal	Polarity:	Positive
Scan type:	Full	Data type:	Centroid

3. Review of chromatography

- Are the hydrophilic peptides retaining as well as the hydrophobic peptides?
- What is the peak width and symmetry?
- Are all expected ions present?
- Is the intensity appropriate to amount injected?

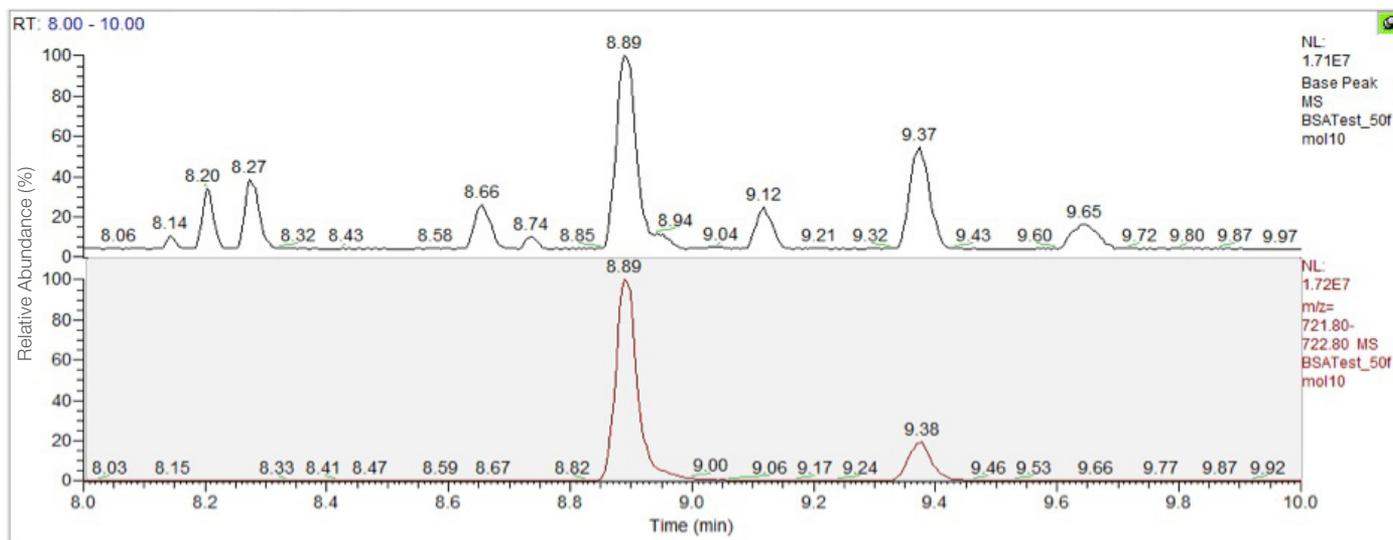
Extraction of known peptide

- Extract a known peptide from the chromatogram, for example *m/z* 722.3 in BSA Digest (YIC*DNQDTISSK)
- Evaluate the peak width and symmetry

Specification for *m/z* 722.3:

PW_{HH} (sec): 1.5–3.5

Asymmetry (EP): 0.9–1.7



4. Post-run cleaning gradients

- After the gradient, add three fast “seesaw” cycles of organic to aqueous
- Review the chromatography – If impurities can be seen eluting during the high organic areas of the “seesaw” gradient, the column is contaminated
- Seesaw gradients can be used to clean the column

Related Thermo Scientific products

88320 Thermo Scientific™ Pierce™ Peptide Retention Time Calibration Mixture

88341 Thermo Scientific™ Pierce™ BSA Protein Digest, MS grade

161089 Thermo Scientific™ Dionex™ Cytochrome C Digest

Current versions of product instructions are available at separatedbyexperience.com/chromexpert.

Learn more about EASY-Spray columns at
thermofisher.com/EASYspray

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representatives for details.

AB21960-EN 0919S

ThermoFisher
SCIENTIFIC