

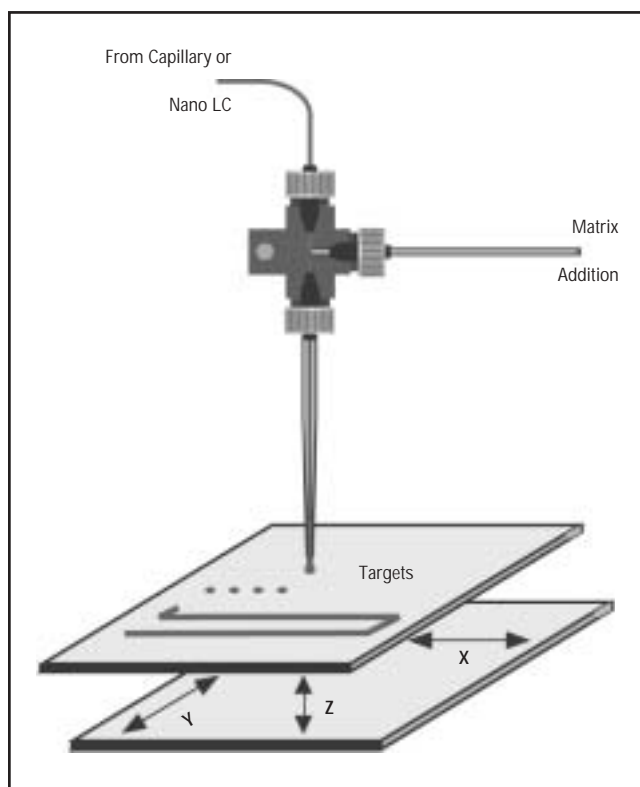
# LC-MALDI MS Using Monolithic Capillary Columns

## INTRODUCTION

Monolithic capillary columns, based on a polystyrene-divinylbenzene polymer (PS-DVB), show excellent separation performance for both proteins and peptides. Using short 5-cm length columns results in very fast protein and peptide separations with a peak width at half height (PWHH) of only a few seconds. Another advantage of the monolithic structure is the very robust column bed, resulting in zero voiding and a superior column lifetime.

## EXPERIMENTAL

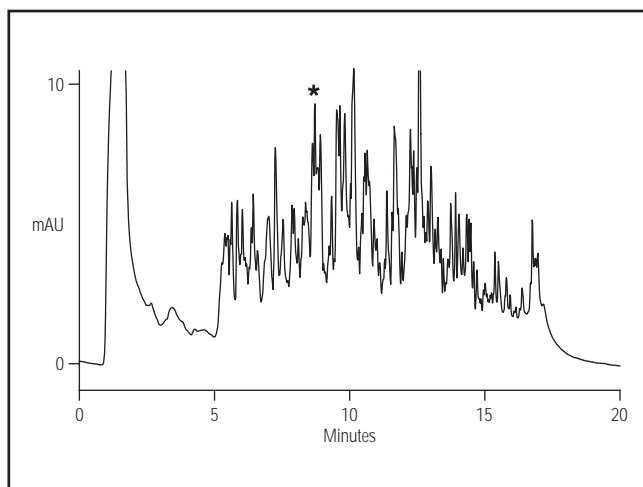
A complex protein digest sample was separated on 200- $\mu\text{m}$  i.d. by a 5-cm Monolithic capillary column using an UltiMate™ Plus Capillary LC System. This capillary LC system was equipped with a capillary flow sensor to assure precise flow delivery and was on-line coupled to a Probot™ Microfraction Collector for subsequent fraction collection. Fractions were collected every 10 s on MALDI targets. Probot's robotic system moves only the collection table and not the collection needle (Figure 1), thus unique precision of  $\pm 2\ \mu\text{m}$  is achieved routinely. During fractionation, an alpha-cyano-4-hydroxy cinnamic acid (CHCA) solution was added 1:1 to the effluent as matrix solution. The fractionated peptides were further analyzed by MALDI-TOF MS.



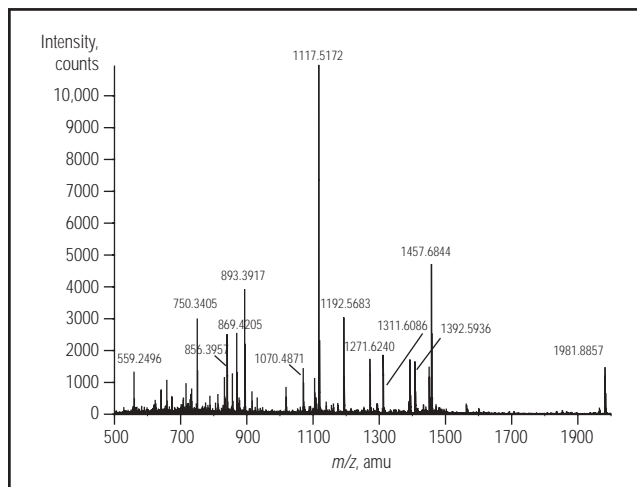
**Figure 1.** X, Y, Z table movement with static needle for high-precision sample spotting on MALDI targets.

Now sold under the  
Thermo Scientific brand

**Thermo**  
SCIENTIFIC



**Figure 2.** UV trace of a digested protein mixture separated on a 200- $\mu\text{m}$  i.d.  $\times$  5 cm Monolithic capillary column and subsequent on-line fraction collection on MALDI targets.



**Figure 3.** MS data from spot #41 (marked with an asterisk in Figure 2).

Figure 2 shows the separation of the digested proteins. A gradient from 0 to 70% acetonitrile is performed in 10 min, resulting in a rapid separation of the peptides. Individual peaks were collected on a high-density MALDI target for subsequent MS analysis. Figure 3 shows the MS spectrum of an arbitrarily chosen peak marked with an asterisk (spot #41).

Coupling Monolithic capillary columns to MALDI-MS results in very fast and sensitive LC-MS analysis, making these columns ideally suited for high-throughput LC-MS proteomics.

### CONDITIONS

LC system:	UltiMate Plus (P/N 163628) FAMOS™ (P/N 163654) Probot (P/N 160560)
Loop:	1 $\mu\text{L}$ (full-loop injections)
Column:	Monolithic capillary column (P/N 161409) 200- $\mu\text{m}$ i.d. $\times$ 5 cm
Flow rate:	2 $\mu\text{L}/\text{min}$
UV-Vis wavelength:	214 nm
Mobile phase:	(A) $\text{H}_2\text{O}$ – 0% acetonitrile + 0.05% TFA; (B) $\text{H}_2\text{O}$ – 50% acetonitrile + 0.04% TFA
Gradient:	0–70% B in 10 min
Matrix solution:	10 mg/mL alpha-cyano-4-hydroxy cinnamic acid (CHCA)
Sample	Tryptic digest of 17 proteins



FAMOS, Probot, and UltiMate are trademarks of Dionex Corporation.

**Dionex Corporation**  
1228 Titan Way  
P.O. Box 3603  
Sunnyvale, CA  
94088-3603  
(408) 737-0700

**Dionex Corporation**  
Salt Lake City Technical Center  
1515 West 2200 South, Suite A  
Salt Lake City, UT  
84119-1484  
(801) 972-9292

**Dionex U.S. Regional Offices**  
Sunnyvale, CA (408) 737-8522  
Westmont, IL (630) 789-3660  
Houston, TX (281) 847-5652  
Atlanta, GA (770) 432-8100  
Marlton, NJ (856) 596-0609

**Dionex International Subsidiaries**  
Australia 61 (2) 9420 5233 Austria (01) 616 51 25 Belgium (03) 353 42 94 Canada (905) 844-9650 China (852) 2428 3282  
Denmark 36 36 90 90 France 01 39 30 01 10 Germany 06126-991-0 Italy (06) 66 51 50 52 Japan (06) 6885-1213  
Korea 82 2 2653 2580 The Netherlands (0161) 43 43 03 Switzerland (062) 205 99 66 United Kingdom (01276) 691722  
\* Designed, developed, and manufactured under an NSAI registered ISO 9001 Quality System.

