

1599-9835

Author:
PerkinElmer, Inc.
Waltham, MA

A New, Sensitive Method for Lignan Metabolite Detection of Flaxseed- Fed Mice Using PerkinElmer QSight® 220

Introduction

Dietary phytochemicals such as polyphenols, carotenoids, saponins and glucosinolates have been suggested to play important roles in human health and may contribute to the prevention of diseases. While the intrinsic food

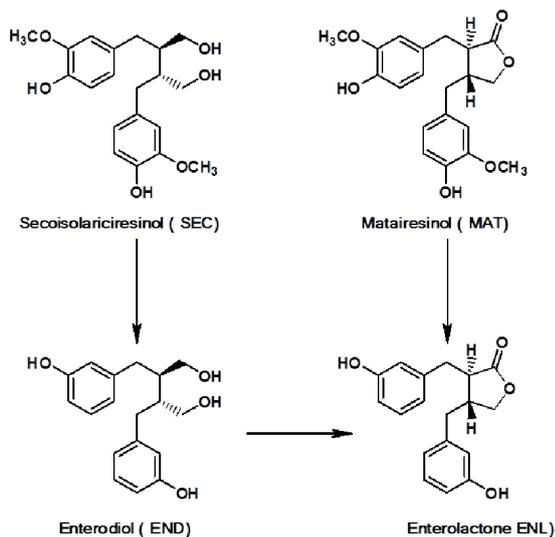
bioactives have been found to act as strong antioxidants and exert various activities *in vitro*, the actual bioactive forms *in vivo* are not clear. In fact, it is often the metabolites of a phytochemical that exert bioactivities. Among the various bioactive polyphenols found in plant-based foods, lignans are one of the most important subgroups. For example, plant lignans such as secoisolarisresinol diglucoside (SDG) glycoside in flax seed are not free molecules but part of the lignan complex. These lignan glycosides are released by digestive enzymes and further hydrolyzed to produce the aglycones secoisolariciresinol (SEC) and matairesinol (MAT), which are converted by the microorganisms in human colon to the bioactive mammalian lignans enterodiol (END) and enterolactone (ENL) (Figure 1)^[1]. However, due to low bioavailability, metabolites of lignans are at minute concentrations in biological fluid samples. This presents a significant challenge for food and nutritional scientists using conventional detectors coupled with liquid chromatography. LC-MS/MS proves beneficial.

2. Method

2.1. Sample Clean Up

All mouse serum samples were taken from a feeding trial with a diet containing different amounts and fractions of flaxseed content. The serum (50 mL) was mixed with 100 mL freshly prepared hydrolysis reagent (contained 2 mg/mL b-glucuronidase/sulfatase in 50 mM NaOAc buffer, pH 4) and incubated at 37 °C for 19 h. 600 mL MeOH was added to each sample (80% final MeOH concentration) after hydrolysis. The samples were vigorously vortexed and placed on a shaker for 1 h, then centrifuged at 20,000 g for 10 min at 4 °C. The supernatant was transferred to a new microfuge tube, and the pellet was washed with 400 mL of 80% MeOH and then centrifuged. The supernatant was combined and evaporated to dryness using the Speed-Vac with solvent compatible ultra-cold trap and vacuum pump. The dried samples were re-dissolved into 50 mL of 80% MeOH by vortexing, sonication for 10 min and centrifuged at 20,000 g for 10 min at 4°C. The supernatant was transferred into insert of HPLC vials and analyzed using the QSight® 220 triple quadrupole.

Figure 1: Flax Lignans & Mammalian Metabolites



2.2. Recovery of Lignans

To verify the accuracy of the sample preparation procedure, recovering was determined by calculating the percent recovery of known amounts of lignan standards added to serum from mouse which was on basal diet. Serum was spiked with 10 mL of 5 mg/mL of lignan standard mixture (MAT, SEC, END and ENL) and hydrolyzed and extracted the same way as unspiked samples

Quick Facts:

- High sensitivity method for quantitation of dietary polyphenols and their metabolites in biological systems using the QSight® 220 triple quadrupole mass spec.
- Analysis of Secoisolariciresinol, Enterodiol, Matairesinol and Enterolactone.
- $R^2 > 0.993$ for all calibration curves.
- Detection limit < 1 ppb for all investigated dietary polyphenols and their metabolites.

2.3. Mass Spectrometry Conditions

The LC-MS/MS analysis was performed using the QSight® 220 triple quadrupole mass spectrometer. Table 1 outlines the parameter settings used during this method.

Table 1: Overview of Settings used on the QSight® 220 Mass Spectrometer

ESI Voltage (V)	5800
HSID Temp (°C)	250
Nebulizer Gas Setting	300
Drying Gas Setting	125
Heating Gas	350
Source Temp. (°C)	250
Dwell Time (ms)	100
Pause Time (ms)	5

2.4. LC Conditions

This method utilized a Shimadzu® Prominence UFLC® XR system. Sample injections of 5 µL were loaded onto an Phenomenex Kinetex® Phenyl-Hexyl 10A column (100x4.6mmx2.6µm) with a KrudKatcher ultra in-line filter as a guard column. Using the gradient flow as shown below in Table 2 at a flow rate of 0.4 mL/min. The mobile phase consisted of solvent A (0.1% formic acid + 99.9% H₂O) and solvent B (0.1% formic acid + 99.9% MeOH).

Table 2: Optimized MRM Parameters

Compound name	Precursor	Fragment	CCL2	CE
Secoisolariciresinol	360.66	164.76	46	32
Enterodiol	300.67	252.78	46	26
Matairesinol	356.62	82.87	40	32
Enterolactone	296.68	252.80	46	26

Table 3: LC Cycle Time

Time (min)	Procedure Solvent B composition (%)
0-10	58
12	80
13	100
16	58
23	58

3. Results

3.1. Lignan Recovery

The conventional method for extraction and cleanup for serum samples include enzyme hydrolysis, liquid/liquid partition, evaporation and solid phase extraction chromatography. Those procedures are often time consuming and costly. This method uses microcentrifugation with Speed-vac and proved to be a simple and valid effective protocol with good recovery rates for the compounds of interest. Table 4 shows the recovery rates for the compounds of interest, proving the method protocol to be effective.

Table 4: Recovery Rate by Standard Addition

Lignan	END	ENL	MAT	SEC
Before Std Addition (ng/mL)	0.7 ± 0.4	3.5 ± 2.4	0.0 ± 0.0	0.6 ± 0.6
After Std Addition (ng/mL)	549.1 ± 36.8	500.8 ± 16.8	528.10 ± 46.90	606.7 ± 22.0
Difference (ng/mL)	548.4 ± 37.1	497.3 ± 19.1	528.1 ± 46.9	606.1 ± 21.9
Expected Difference (ng/mL)	500.0	500.0	500.0	500.0
Recovery Rate %	109.7 ± 7.4	99.5 ± 3.8	105.6 ± 9.4	121.2 ± 4.4

The recovery rates for these compounds were between 99-121% as determined by standard addition of known amounts of serum from mouse fed basal diet.

3.2. Detection of Lignans at Low Concentration

All the lignans and their metabolites were well-separated and detected using the QSight® 220 triple quadrupole mass spectrometer. Figure 2 & 3 illustrates the results for a typical LC-MS/MS total ion chromatogram of the standards (5ppb=15nM) and actual serum sample respectively. Identification of the lignans was done by congruent retention time and (-)-ESI mass spectrometry. For sensitive quantification, the MS/MS pairs were between the molecular ion [M-H]⁻¹ and its most abundant daughter ion.

Figure 2: Example Total Ion Chromatogram of the Standards at 5ppb=15nM

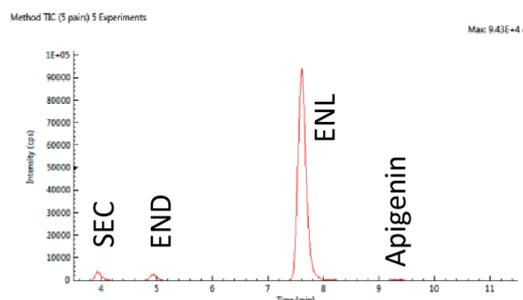
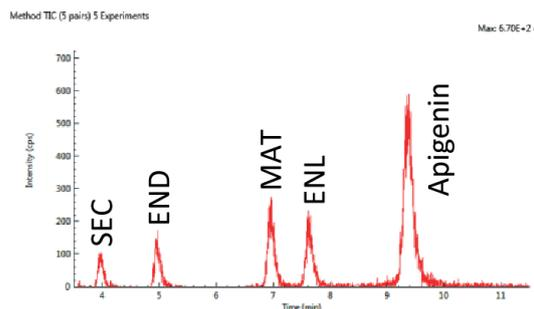


Figure 3: Example Total Ion Chromatogram of the Actual Serum Sample.



Using this method, the detection of all targeted lignans (SEC and MAT) and the two most important metabolites (END and ENL) in the serum samples of mice fed different amount of flax seed meals was detected. ENL and END are mammalian lignans known for their many health beneficial effects. In addition, a minute amount of apigenin, a flavonoid also found in flaxseed was detected in the serum (Figure 2 & 3).

Table 5: Bioavailability of Lignans & Metabolites in Healthy Male Mice Fed Different Flax Seed Fractions (μM in serum) (selected samples).

Treatment	sample	END μM	ENL μM	MAT μM	SEC μM
Basal diet (control)	S1	0.000	0.007	0.000	0.000
	S2	0.000	0.006	0.000	0.000
10% Flaxseed	S7	4.324	3.541	0.000	0.298
	S8	2.983	0.678	0.000	0.088
6% Flax kernel	S9	0.071	0.058	0.000	0.070
	S10	0.049	0.088	0.000	0.000
4% Flax hull	S31	3.374	1.678	0.000	0.082
	S32	2.938	0.247	0.000	0.023

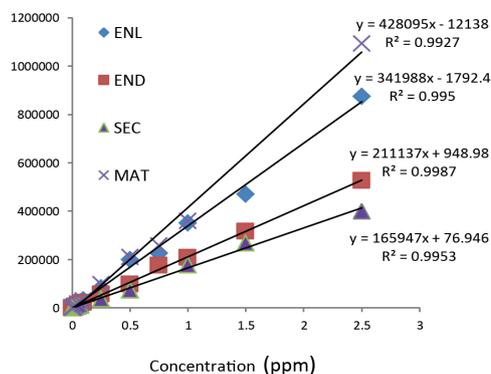
Table 5 shows the serum concentrations of the four lignans in mice fed diets containing no or different flaxseed components, i.e. 10% whole flaxseed flour, 6% flax kernel flour or 4% flax hull powder. All these compounds were detected at the nM – μM level. There were individual differences in the ability to produce the metabolites, possibly due to the different microbial flora of these mice. Some produced more END than ENL, others more ENL than END. The two plant lignan aglycones SEC and MAT were of very low bioavailability, in fact, no MAT was detected in the majority of the samples. These plant lignans are mostly in the hull of the flax seed, therefore, it is highly significant to see the low concentrations of END and ENL in serum of mice fed kernel.

The concentrations of END, ENL, SEC and MAT were measured for 180 samples and they differed significantly, with END to be the most dominant metabolite ranging from 0.00452 to 100.35 μM , followed by ENL (0.00452 to 61.75 μM), SEC (0.0048 to 49.45 μM), and MAT (0.0014 to 2.76 μM). Enterodiol (END), a mammalian lignan, was the predominant metabolite in the serum whereas the plant lignan matairesinol (MAT), was found to be barely absorbed.

3.3. Linearity

The standard curves around the concentrations detected in the serum samples were highly linear for all analytes, with $R^2 > 0.993$ with no weighting (Figure 4). The detection limits for all compounds were $< 1 \text{ ppb}$, ca. 3 nM.

Figure 4: Standard Curves of Flax Lignans and their Metabolites



The calibration curves generated for SEC (360.66/164.76), END (300.67/252.78), MAT (356.62/82.87) and ENL (296.68/252.80) showed good linearity ($R^2 > 0.993$).

4. Conclusion

Present results demonstrate that the reported LC-MS/MS method using a new highly sensitive PerkinElmer QSiht[®] 220 triple quadrupole platform is capable of detecting $< 1 \text{ ppb}$ level of dietary polyphenols and their metabolites in biological systems using only 50 μL of biological sample. This new and robust analytical method can help better understand the health benefits and risks of lignan containing foods.

5. References

- Sok, DE; Cui, HS; Kim, MR. Isolation and Bioactivities of Furfuran Type Lignan Compounds from Edible Plants. Recent Patents on Food, Nutrition & Agriculture, 2009, 1, 87-95
- Wu, Q; Wang, M; Simon, JE. Analytical methods to determine phytoestrogenic compounds. Journal of Chromatography B, 812 (2004) 325–355

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

PerkinElmer, Inc.
Wallac Oy
PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357

ISO 13485
ISO 9001
CMDCAS
ISO 14001
OHSAS 18001



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2018, PerkinElmer, Inc. All rights reserved. PerkinElmer[®] is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

1599-9835