

LC-MS/MS 法测定辛伐他汀的药物一致性评价方法学考察

王立杰 蔡悠悠

赛默飞世尔科技（中国）有限公司

摘要

建立一种测定人体血浆中辛伐他汀的 LC-MS/MS 方法，辛伐他汀 d6 作为内标，采用乙腈沉淀蛋白，离心后上清液用水复溶的前处理方法。分析柱为 Thermo Hypersil Gold (100 mm*2.1 mm, 3 μ) 色谱柱，用 PH 值为 4.0 的 2.00 mM 醋酸铵水溶液与乙腈梯度洗脱。离子源为可加热电喷雾离子源（HESI），正离子检测，扫描方式为多反应监测（SRM），辛伐他汀和内标的监测离子对分别为 m/z 436.2 $\rightarrow$ 199.1 和 442.3 $\rightarrow$ 199.1。分析时间为 5.5 min，辛伐他汀在 0.05-50 ng/mL 浓度范围内线性良好。辛伐他汀定量下限为 0.05 ng/mL。本方法的灵敏度、精确度和准确度、基质效应、回收率等符合生物样品分析要求，本方法符合法规要求，可以用于药物生物等效性一致性评价的研究。

关键词

辛伐他汀；LC-MS/MS；一致性评价

1. 引言

辛伐他汀是由土曲霉内提炼而得，最早是由默克药厂所开发，并于 1992 年开始进入医疗用途。该药列名于世界卫生组织基本药物清单，属于基础医疗体系必备药物之一^[1]。作为一种口服降血脂药物，常见商品名“Zocor”。该品会在运动、节食，和减肥时服用，以避免发生血脂升高。此外辛伐他汀也可降低高心脏病风险者发作的机会。严重副作用包含肌病变、肝脏伤害，以及血糖升高。常见副作用则有便秘、头痛，和恶心。有肾脏问题者宜降低剂量。有证据显示妊娠期间服用该药物会对胎儿造成伤害，哺乳期间不得使用该品。辛伐他汀属于 HMG-CoA 还原酶抑制剂，借此抑制肝脏制造胆固醇。鉴于辛伐他汀的广泛用途，越来越多的生产厂开始生产或仿制辛伐他汀制剂。为了方便进行辛伐他汀的生物等效性一致性评价，本文建立一种灵敏、快速、可靠的 LC-MS/MS 检测方法，并按中国药典按照（CFDA）的指导原则进行了方法学验证。

2. 实验方法

2.1 试剂与样品制备

辛伐他汀（纯度 98%）购自 Toronto Research Chemicals，HPLC 级甲醇、乙腈购自 Fisher Chemical，超纯水自制，其他试剂为分析级。

2.2 实验用人血浆

购买自 Bioreclamation LLC。

2.3 实验仪器

分析仪器为液相色谱串联三重四极杆质谱系统，其中液相色谱为 Vanquish Binary Horizon HPLC，质谱为 TSQ Quantis（ThermoFisher Scientific），三重四极杆配有可加热电喷雾离子源。Thermo Hypersil Gold (100mm*2.1mm,3 μ) 为分析柱。

2.4 LC-MS/MS 条件

液相条件：流动相为 PH 值为 4.0 的 2.00 mM 醋酸铵水溶液和乙腈，梯度洗脱，流速为 0.4 ml/min，分析时间为 5.5 min。

MS/MS 条件：喷雾电压为 4000V，离子源和离子传输管温度分别为 350℃和 350℃；碰撞气氦气压力为 1.5 mTorr；辛伐他汀和内标的监测离子对分别为 m/z 436.2 → 199.1 和 442.3 → 199.1。

2.5 储备液配置和标准曲线建立

辛伐他汀的储备液浓度为 0.5 mg/mL，溶剂为乙腈。将辛伐他汀储备液稀释成浓度为 12500, 2500, 125, 100, 50, 25, 5 ng/mL 的工作溶液。

2.6 血浆处理

100mL 空白血浆加入 100mL 200mM 乙酸铵水溶液，用乙酸调 PH 至 4.5，备用。

处理后的空白血浆中加入辛伐他汀标准溶液和内标溶液，制备高、中、低质控样品。

2.7 方法学验证

按照 CFDA 的指导原则进行方法学验证。

2.7.1 专属性

通过对比空白血浆、加入内标溶液的空白血浆和加入辛伐他汀配制低质控浓度的血浆样品谱图，计算低质控样品的准确度，考察该方法的专属性。

2.7.2 最低定量限

以标准曲线最低浓度点作为质控样品，计算其准确度考察最低定量限。

2.7.3 建立标准曲线

取 95 μL 空白血浆，加入 5 μL 辛伐他汀标准工作溶液和内标溶液，经样品前处理后进样分析。分别以辛伐他汀的峰面积与内标峰面积的比值对其浓度作图，用加权最小二乘法进行线性回归，得加权标准曲线。

2.7.4 精密度和准确度

高、中、低质控浓度 6 个样品一日内进样，计算日内精密度和准确度。每个样品连续 3 日内分别进样分析，计算日间精密度和准确度。

2.7.5 回收率和基质效应

血浆样品在沉淀蛋白前和沉淀蛋白后分别加入辛伐他汀标准溶液和内标溶液，计算沉淀前后的比值考察本方法的回收率。在血浆沉淀蛋白后的上清液和 60% 乙腈水溶液中分别加入辛伐他汀标准溶液和内标溶液，计算两者的比值考察本方法的基质效应。

2.7.6 稳定性实验

辛伐他汀储备液、内标储备液、最高与最低两个浓度工作溶液经 4℃保存 15 天，计算 4℃保存的溶液与新鲜配制的溶液之间的偏差考察溶液的长期稳定性。

高、低两个质控浓度样品经反复冻融、湿冰保存 2h 和 -80℃保存 15 天、45 天，计算质控样品的准确度考察短期与长期稳定性。

经过前处理后的同一序列间隔 96h 分别进样，考察自动进样器的稳定性。

2.7.7 稀释效应

将 250 ng/mL 辛伐他汀样品用空白血浆稀释 10 倍。计算准确度考察稀释对样品测定的影响。

2.8 高脂血效应

用高脂血浆配制高、低质控样品，计算准确度考察高脂血效应。

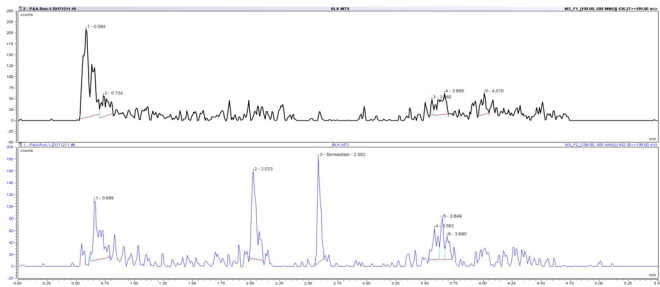
2.9 序列大小

计算标准曲线与质控样品的准确度，考察一个序列中最大容纳的样品数量。

3. 实验结果与讨论

3.1 专属性

通过对比空白血浆、加入内标溶液的空白血浆和加入辛伐他汀配制低质控浓度的血浆样品色谱图，并低质控浓度样品的准确度，结果如下。说明血浆中内源性物质对辛伐他汀和内标的检测没有干扰。



Calibration Curve Parameters for Simvastatin in Human plasma

Analysis Group	Slope	y-Intercept	Correlation Coefficient
001	55.64169183	-0.08225152632	0.9956
002	55.01285927	0.294891413	0.9974
003	53.47304402	0.833452586	0.9954

Calibration Curve Data for Simvastatin in Human plasma

Theoretical Concentration (ng/mL)								
Analysis Group	0.05	0.10	0.50	2.0	10.0	25.0	45.0	50.0
001	0.053	*0.120	0.450	2.017	9.469	24.606	44.949	50.751
	0.044	0.111	0.480	2.117	10.009	24.802	45.415	52.846
002	0.050	*0.083	0.472	1.872	10.370	24.788	46.422	49.323
	0.053	0.089	0.520	1.970	10.407	25.146	45.394	52.321
003	0.046	0.118	0.501	2.104	9.891	24.573	45.424	51.926
	*0.034	0.100	0.499	1.869	9.440	25.251	45.054	47.543
n	5	4	6	6	6	6	6	6
Mean	0.0492	0.106	0.485	2.02	10	24.8	45.5	51.4
SD	0.00409	0.0151	0.0269	0.101	0.385	0.228	0.542	1.41
RSD (%)	8.3	14.2	5.5	5.0	3.9	0.9	1.2	2.7
Accuracy (%)	98.4	106.0	97.0	101.0	100.0	99.2	101.1	102.8

* Failed acceptance criteria

3.3 最低定量限

辛伐他汀样品的最低定量限为 0.05 ng/mL。

Sensitivity Data for Simvastatin in Human plasma

Analysis Group	Theoretical Concentration (ng/mL)	Calculated Concentration (ng/mL)	Diff (%)
001	0.05	0.052	4.0
		0.057	14.0
		0.061	22.0
		0.053	6.0
		0.053	6.0
		0.058	16.0
n		6	
Intra-Assay Mean		0.0557	
SD		0.00356	
RSD (%)		6.4	
Accuracy (%)		111.4	

002	0.05		0.037	-26.0
			0.054	8.0
			0.053	6.0
			0.042	-16.0
			0.05	0.0
			0.049	-2.0
n		6		
Intra-Assay Mean		0.0475		
SD		0.00666		
RSD (%)		14.0		
Accuracy (%)		95.0		
00.3	0.05		0.042	-16.0
			0.034	-32.0
			0.055	10.0
			0.057	14.0
			0.056	12.0
			0.052	4.0
n		6		
Intra-Assay Mean		0.0493		
SD		0.00929		
RSD (%)		18.8		
Accuracy (%)		98.6		
Overall				
n		18		
Inter-Assay Mean		0.0508		
SD		0.00742		
RSD (%)		14.6		
Accuracy (%)		101.6		

3.4 精密度和准确度

高、中、低浓度质控样品的日间和日内精密度和准确度均小于 15%，低浓度小于 20%，符合生物样品分析要求。

Quality Control Sample Data for Simvastatin in Human plasma

Analysis Group		Theoretical Concentration (ng/mL)					
001		0.15	Diff (%)	2.50	Diff (%)	37.50	Diff (%)
		0.168	12.0	2.542	1.7	41.650	11.1
		0.184	22.7	2.408	-3.7	43.271	15.4
		0.123	-18.0	2.545	1.8	39.711	5.9
		0.167	11.3	2.659	6.4	41.425	10.5
		0.164	9.3	2.698	7.9	42.934	14.5
		0.169	12.7	2.656	6.2	42.339	12.9
	n	6		6		6	

Intra-Assay Mean	0.163		2.58		41.9	
SD	0.0206		0.108		1.28	
RSD (%)	12.6		4.2		3.1	
Accuracy (%)	108.7		103.2		111.7	
002	0.146	-2.7	2.744	9.8	42.705	13.9
	0.148	-1.3	2.752	10.1	43.630	16.3
	0.198	32.0	2.716	8.6	41.227	9.9
	0.146	-2.7	2.804	12.2	43.115	15.0
	0.149	-0.7	2.791	11.6	40.796	8.8
	0.169	12.7	2.735	9.4	44.287	18.1
n	6		6		6	
Intra-Assay Mean	0.159		2.76		42.6	
SD	0.0209		0.0338		1.37	
RSD (%)	13.1		1.2		3.2	
Accuracy (%)	106.0		110.4		113.6	
003	0.145	-3.3	2.603	4.1	37.779	0.7
	0.164	9.3	2.627	5.1	39.645	5.7
	0.140	-6.7	2.718	8.7	39.1	4.3
	0.171	14.0	2.800	12.0	39.62	5.7
	0.166	10.7	2.759	10.4	39.212	4.6
	0.164	9.3	2.856	14.2	40.067	6.8
n	6		6		6	
Intra-Assay Mean	0.158		2.73		39.2	
SD	0.0126		0.0984		0.793	
RSD (%)	8.0		3.6		2.0	
Accuracy (%)	105.3		109.2		104.5	
Overall						
n	18		18		18	
Inter-Assay Mean	0.16		2.69		41.3	
SD	0.0174		0.112		1.86	
RSD (%)	10.9		4.2		4.5	
Accuracy (%)	106.7		107.6		110.1	

3.5 回收率和基质效应

辛伐他汀低、中、高质控样品的回收率分别在 98.2%、96% 和 91.1%，内标低、中、高质控样品的回收率为 90.8%、88.1% 和 87.3%，说明本实验前处理方法合适。辛伐他汀低、高质控样品经过内标归一化的基质效应分别为 0.981 和 1.000，说明本方法没有显著的基质效应。

Recovery Data for Simvastatin and the Internal Standard in Human plasma

Theoretical Concentration (ng/mL)	Simvastatin		Internal Standard	
	Extracted QC Sample Peak Area	Recovery Sample Peak Area	Extracted QC Sample Peak Area	Recovery Sample Peak Area
0.15	999	945	12029	12579
	1028	1115	12205	13100
	1197	1136	10692	12663
	944		11339	
	1007		11887	
	1104		11482	
n	6	3	6	3
Mean	1046.5	1065.3	11605.7	12780.7
SD	90.1	104.7	554.8	279.7
RSD (%)	8.6	9.8	4.8	2.2
Recovery (%)		98.2		90.8
2.5	17637	17226	11660	12670
	17319	18370	11419	12885
	17088	18739	11415	13436
	16727		10823	
	17884		11624	
	17708		11745	
n	6	3	6	3
Mean	17393.8	18111.7	11447.7	12997.0
SD	434.0	788.9	333.7	395.1
RSD (%)	2.5	4.4	2.9	3.0
Recovery (%)		96.0		88.1
37.5	266419	306422	11339	13796
	278968	299496	11621	13038
	274490	304232	12101	13685
	281748		11877	
	272109		12123	
	285356		11711	
n	6	3	6	3
Mean	276515.0	303383.3	11795.3	13506.3
SD	6884.9	3540.1	301.1	409.4
RSD (%)	2.5	1.2	2.6	3.0
Recovery (%)		91.1		87.3
Overall Recovery (%)		95.1		88.7

Matrix Factor Data for Simvastatin and the Internal Standard in Human plasma

Simvastatin			Internal Standard						
Matrix ID	Post Extraction Spike Peak Area	Pure Solution Peak Area	Mean Pure Solution Peak Area	Matrix Factor	Post Extraction Spike Peak Area	Pure Solution Peak Area	Mean Pure Solution Peak Area	Matrix Factor	ISTD Normalized Matrix Factor
Lot1	918.0	1090.0	1111.7	0.826	9930.0	13171.0	13052.7	0.761	1.09
Lot2	789.0	1103.0		0.710	9594.0	12972.0		0.735	0.966
Lot3	756.0	1142.0		0.680	9164.0	13015.0		0.702	0.969
Lot4	827.0			0.744	9502.0			0.728	1.02
Lot5	745.0			0.670	9598.0			0.735	0.912
Lot6	754.0			0.678	9545.0			0.731	0.927
n				6				6	6
Mean				0.718				0.732	0.981
SD				0.0596				0.0188	0.0655
RSD (%)				8.3				2.6	6.7
Lot1	217534.0	304163.0	293171.7	0.742	9578.0	12829.0	12389.0	0.734	1.01
Lot2	218332.0	289525.0		0.745	9133.0	12318.0		0.700	1.06
Lot3	219329.0	285827.0		0.748	9731.0	12020.0		0.746	1.00
Lot4	225575.0			0.769	10225.0			0.783	0.982
Lot5	217642.0			0.742	10051.0			0.770	0.964
Lot6	213755.0			0.729	9500.0			0.728	1.00
n				6				6	6
Mean				0.746				0.744	1.00
SD				0.0131				0.03	0.0325
RSD (%)				1.8				4.0	3.3

3.6 稳定性

辛伐他汀储备液、内标储备液、最高与最低两个浓度工作溶液在 4℃ 保存 15 天的条件下稳定

Solution Stability for Simvastatin Stored at 4℃ for 15 Days

Time	Stock Standard Solution		Internal Standard		Lowest Concentration Working Solution		Highest Concentration Working Solution	
	0 Days	15 Days	0 Days	15 Days	0 Days	15 Days	0 Days	15 Days
Peak Area Ratio	2.4779	2.3844	0.4036	0.4194	0.0295	0.0271	1.2133	1.1432
Peak Area Ratio	2.4440	2.3874	0.4092	0.4189	0.0293	0.0270	1.2139	1.1494
Peak Area Ratio	2.4665	2.4006	0.4054	0.4166	0.0293	0.0267	1.2178	1.1496
Peak Area Ratio	2.4706	2.4095	0.4048	0.4150	0.0293	0.0264	1.2161	1.1514
Peak Area Ratio	2.4800	2.4077	0.4032	0.4153	0.0296	0.0269	1.2113	1.1506
Peak Area Ratio	2.4865	2.4093	0.4022	0.4151	0.0290	0.0272	1.2157	1.1427
n	6	6	6	6	6	6	6	6
Mean PeakArea Ratio	2.47	2.40	0.405	0.417	0.0293	0.0269	1.21	1.15
Combined Mean	2.44		2.96		-8.2		-4.96	

Short-Term Matrix Stability Data for Simvastatin in Human plasma
After Storage at -80℃ for 15 Days

Theoretical Concentration (ng/mL)	0.15	Diff (%)	37.5	Diff (%)
	0.156	4.0	40.358	7.6
	0.161	7.3	40.952	9.2
	0.176	17.3	40.714	8.6
	0.149	-0.7	35.808	-4.5
	0.160	6.7	37.420	-0.2
	0.173	15.3	38.238	2.0
n	6		6	
Mean	0.163		38.9	
SD	0.0103		2.09	
RSD (%)	6.3		5.4	
Accuracy (%)	108.7		103.7	

After Storage at -80℃ for 45 Days

Theoretical Concentration (ng/mL)	0.15	Diff (%)	37.5	Diff (%)
	0.154	2.7	31.011	-17.3
	0.164	9.3	32.960	-12.1
	0.164	9.3	34.333	-8.4
	0.142	-5.3	34.346	-8.4
	0.129	-14.0	30.482	-18.7
	0.180	20.0	35.328	-5.8
n	6		6	
Mean	0.156		33.1	
SD	0.0181		1.96	
RSD (%)	11.6		5.9	
Accuracy (%)	104.0		88.3	

经过前处理后的样品序列在自动进样器中存放 96h 稳定性良好。

Reinjection Reproducibility Data for Simvastatin in Human plasma
Storage at 4℃ for 96 Hours

Theoretical Concentration (ng/mL)	Calculated Concentration (ng/mL)		Diff (%)
0.15	0.156		4.0
	0.165		10.0
	0.166		10.7
	0.167		11.3
	0.155		3.3
	0.170		13.3
Mean	0.163		108.7
2.5	2.618		4.7
	2.581		3.2
	2.867		14.7
	2.730		9.2
	2.721		8.8
	2.776		11.0
Mean	2.72		108.8
37.5	41.46		10.6
	39.806		6.1
	40.248		7.3
	39.148		4.4
	41.733		11.3
	40.487		8.0
Mean	40.5	Accuracy (%)	108.0
Original Analysis Group:	3	Reinjection Analysis Date:	96 hours
Slope:	53.47304402	Slope:	53.55975644
y-Intercept:	0.833452586	y-Intercept:	1.007984204
Correlation Coefficient:	0.9954	Correlation Coefficient:	0.9957

3.7 稀释效应

辛伐他汀 250ng/mL 浓度样品经 10 倍体积稀释后测定其准确度和精密度分别为 9.9% 和 3.8%，符合法规要求。

Sample Dilution Analysis for Simvastatin in Human plasma

Theoretical Concentration (ng/mL)	10X Dilution	Diff (%)
250	23.807	-90.5
	24.004	-90.4
	25.778	-89.7
	25.394	-89.8
	24.139	-90.3
	25.858	-89.7
n	6	
Mean	24.8	
SD	0.947	
RSD (%)		3.8
Accuracy (%)	9.9	

3.8 高脂血效应

辛伐他汀低、高浓度高脂血样品测定准确度分别为 102% 和 104.8%，符合法规要求

Assessment for the Quantitation of Simvastatin in Hyperlipidaemic Human plasma

Theoretical Concentration	0.15 (ng/mL)	37.5 (ng/mL)
	0.154	39.789
	0.147	39.196
	0.160	40.651
	0.147	37.758
	0.154	38.068
	0.157	40.528
n	6	6
Mean	0.153	39.3
SD	0.00527	1.22
RSD (%)	3.4	3.1
Accuracy (%)	102.0	104.8

3.9 序列大小

经测定一个序列中最大可容纳 192 个样品。

4. 结论

本文建立了一种灵敏、快速的测定辛伐他汀的 LC-MS/MS 方法，经方法学验证本方法的专属性、精密度和准确度、基质效应、回收率和稳定性等符合生物样品分析法规要求，适用于药物生物等效性一致性评价研究。



赛默飞
官方微信



赛默飞色谱
和质谱中国

热线 800 810 5118
电话 400 650 5118
www.thermofisher.com

ThermoFisher
SCIENTIFIC