



Estimation of Ethylene Glycol and Diethylene Glycol in Propylene Glycol, Glycerin, and Syrup Samples with the Agilent 8890 GC

Author

Praveen Arya and Vivek Dhyani
Agilent Technologies, Inc.

Abstract

This application note demonstrates the use of the Agilent 8890 GC system coupled with an FID detector to detect and quantify ethylene glycol and diethylene glycol in propylene glycol, glycerin, and syrup samples.

The method provides high confidence in results for routine analysis in the pharma industry, whether in production, processing, storage, or commercial testing of syrup samples, or for academic purposes.

Introduction

Propylene glycol and glycerin are commonly used in medicinal syrups as excipients⁵. The solubility of active ingredients is enhanced by their use during formulation⁵. These excipients should be tested for any contamination by ethylene glycol (EG) and diethylene glycol (DEG) as mentioned in regulations such as Indian Pharmacopeia and USP-NF monographs. Some USP-NF monographs include, as part of the applicable identity testing, a limit test for DEG and EG. The relevant safety limit for DEG and EG is not more than (NMT) 0.10%, as recognized by the applicable USP-NF monograph³.

In this application note, we have shared the results of experiments using an Agilent 8890 GC to identify and quantify EG and DEG in matrices like propylene glycol, glycerin, and syrup samples.



Experimental

Instrument parameters

Table 1. Method Parameters

Parameter	Value
Inlet Temperature	250 °C
Inlet Liner	Ultra Inert, low pressure drop, split liner, 4 mm ID (p/n 5190-2295)
Inlet Septa	Inlet septa, long life, 11 mm (p/n 8010-0239)
Injection Volume	0.5 µL
Column	Agilent J&W DB-624 GC column, 30 m × 0.53 mm × 3.00 µm, 7 inch cage, (p/n: 125-1334)
Column Flow	Helium, 2.5 mL/min
Split Ratio	10:1
Oven Program	70 °C for 1 min
	6 °C/min to 150 °C, hold 3 min
	25 °C/min to 245 °C, hold 12 min
FID Temperature	250 °C
FID H2 Flow	40 mL/min
FID Air Flow	300 mL/min
FID Make Up Gas	Nitrogen, 25 mL/min

Standard calibration

Standard solution-1 was prepared by weighing 100 mg each of EG and DEG in a 100 mL volumetric flask and making up the volume to 100 mL with methanol, with thorough mixing. This solution has a concentration of 1,000 µg/mL for EG and DEG.

IS solution-1 was prepared by weighing 100 mg of 1,3-propanediol in a 100 mL volumetric flask and making up the volume to 100 mL with methanol, with thorough mixing. This solution has a concentration of 1,000 µg/mL for 1,3-propanediol.

Using the above standards, the calibration standards were prepared as presented in Table 2.

Table 2. Calibration standards.

	Standard Solution-1, 1,000 µg/mL (mL)	IS Working Solution, 1,000 µg/mL (mL)	Make Up Volume with Methanol (mL)	Final Concentration for EG and DEG (µg/mL)	Final Concentration for 1,3-propanediol (µg/mL)
Calibration Standard-6	5	0.5	10	500	50
Calibration Standard-5	2.5	0.5	10	250	50
Calibration Standard-4	1	0.5	10	100	50
Calibration Standard-3	0.5	0.5	10	50	50
Calibration Standard-2	0.25	0.5	10	25	50
Calibration Standard-1	0.1	0.5	10	10	50

Sample preparation

500 mg of sample was added to a 10 mL volumetric flask. 500 µL of IS solution was added and the volume was made up to 10 mL with methanol, with thorough shaking.

Results and discussion

Standard chromatograms

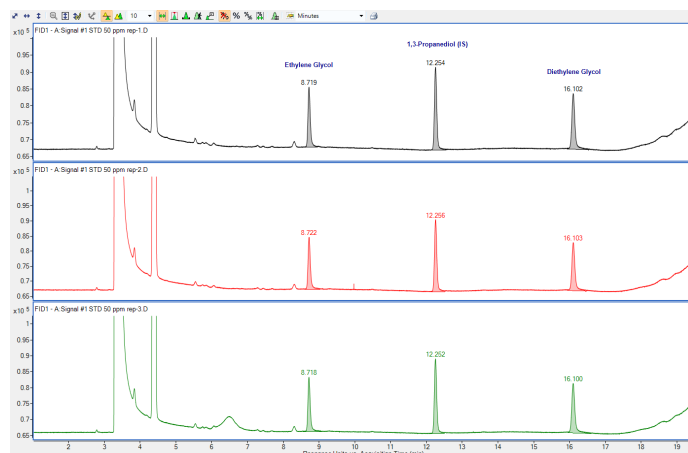


Figure 1. Three replicates of 50 µg/mL EG and DEG standards with IS using methanol as diluent.

Figure 1 shows the overlay of chromatograms obtained from three replicate injections of EG, DEG, and IS (1,3-propanediol) at 50 µg/mL concentration level. The peak shapes, resolution, and repeatability were found to be good. Figures 2 and 3 show the six-point internal-standard-based linearity for calibration standards of EG and DEG, respectively. Good r^2 values were obtained (0.997 for EG and 0.998 for DEG). Table 3 shows the repeatability in terms of internal standard RSD for RT and area for six calibration standards.

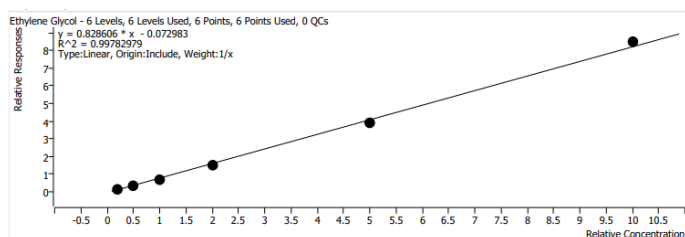


Figure 2. Six-point calibration of EG for 10, 25, 50, 100, 250, and 500 µg/mL, respectively.

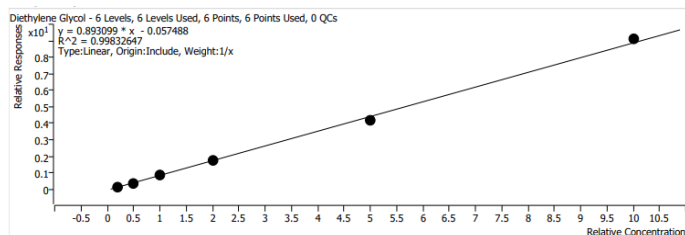


Figure 3. Six-point calibration of DEG for 10, 25, 50, 100, 250, and 500 µg/mL, respectively.

Table 3. %RSD for RT and area for internal standard in six-point calibration standards.

	IS RT (min)	IS Area (Counts)
STD-1_10 µg/mL	12.249	113107
STD-2_25 µg/mL	12.252	109688
STD-3_50 µg/mL	12.255	106848
STD-4_100 µg/mL	12.255	110376
STD-5_250 µg/mL	12.244	117225
STD-6_500 µg/mL	12.255	111914
Mean	12.252	111526.333
SD	0.0045	3513.21
%RSD	0.036	3.15

Table 4. Recovery and repeatability of EG and DEG in 200 µg/mL spiked propylene glycol sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
200 µg/mL spike-1	204.70	102.35	222.05	111.025
200 µg/mL spike-2	203.22	101.61	220.96	110.48
200 µg/mL spike-3	204.86	102.43	214.37	107.185
Mean	204.26	102.13	219.13	109.56
SD	0.904		4.155	
%RSD	0.443		1.896	

Table 5. Recovery and repeatability of EG and DEG in 500 µg/mL spiked propylene glycol sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
500 µg/mL spike-1	456.42	91.284	501.87	100.374
500 µg/mL spike-2	478.86	95.772	502.44	100.488
500 µg/mL spike-3	477.91	95.582	499.81	99.962
Mean	471.06	94.21	501.37	100.27
SD	12.690		1.384	
%RSD	2.694		0.276	

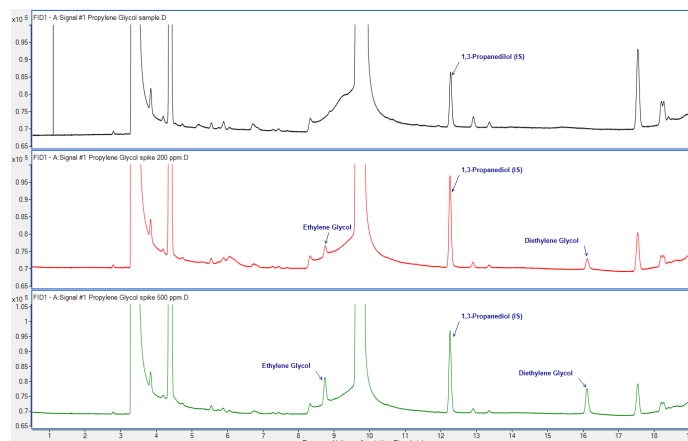


Figure 4. Chromatograms of (from top to bottom) blank propylene glycol sample, and 200 and 500 µg/mL spikes.

Figure 4 shows chromatograms obtained from injection of propylene glycol, and propylene glycol spiked with 200 and 500 µg/mL of EG and DEG, respectively. These chromatograms show good resolution of the compounds of interest from matrix peaks. Tables 4 and 5 show the recovery and repeatability (in terms of %RSD) for propylene glycol sample spiked with 200 and 500 µg/mL of EG and DEG, respectively. Excellent recovery and repeatability were achieved for this experiment with propylene glycol sample.

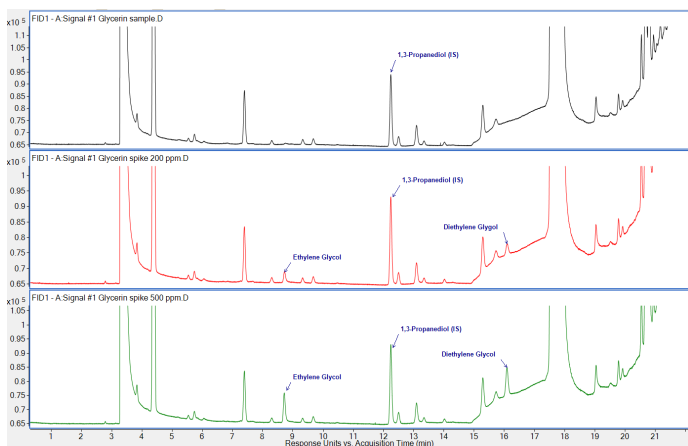


Figure 5. Chromatograms of (from top to bottom) blank glycerin sample, and 200 and 500 µg/mL spikes.

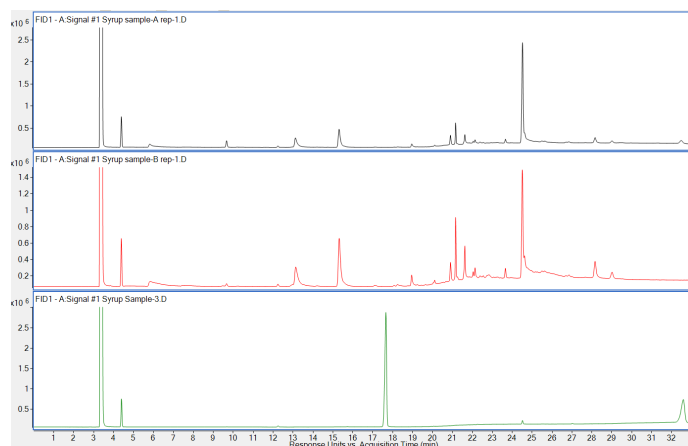


Figure 6. Chromatograms of three different brands of syrup samples.

Table 6. Recovery and repeatability of EG and DEG in 200 µg/mL spiked glycerin sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
200 µg/mL spike-1	243.15	121.575	239.51	119.755
200 µg/mL spike-2	241.32	120.66	244.1	122.05
200 µg/mL spike-3	244.47	122.235	232.55	116.275
Mean	242.98	121.49	238.72	119.36
SD	1.582		5.815	
%RSD	0.651		2.436	

Table 7. Recovery and repeatability of EG and DEG in 500 µg/mL spiked glycerin sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
500 µg/mL spike-1	530.25	106.05	532.81	106.562
500 µg/mL spike-2	536.04	107.208	547.87	109.574
500 µg/mL spike-3	540.16	108.032	547.6	109.52
Mean	535.48	107.10	542.76	108.55
SD	4.978		8.618	
%RSD	0.930		1.588	

Figure 5 shows chromatograms obtained from injection of glycerin, and glycerin spiked with 200 ppm and 500 µg/mL of EG and DEG, respectively. Compounds of interest are well resolved from matrix peaks. Tables 6 and 7 show the recovery and repeatability (in terms of %RSD) for glycerin sample spiked with 200 and 500 µg/mL of EG and DEG, respectively. Excellent recovery and repeatability were achieved for this experiment with glycerin sample.

Table 8. Recovery and repeatability of EG and DEG in 200 µg/mL spiked syrup sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
200 µg/mL spike-1	208.04	104.02	185.17	92.585
200 µg/mL spike-2	214.14	107.07	187.7	93.85
200 µg/mL spike-3	210.9	105.45	181.68	90.84
Mean	211.03	105.51	184.85	92.43
SD	3.052		3.023	
%RSD	1.446		1.635	

Table 9. Recovery and repeatability of EG and DEG in 500 µg/mL spiked syrup sample.

Spiked Sample	EG Results		DEG Results	
	Calculated Amount (µg/mL)	Recovery (%)	Calculated Amount (µg/mL)	Recovery (%)
500 µg/mL spike-1	433.18	86.636	389.61	77.922
500 µg/mL spike-2	429.64	85.928	392.47	78.494
500 µg/mL spike-3	431.82	86.364	385.83	77.166
Mean	431.55	86.31	389.30	77.86
SD	1.786		3.331	
%RSD	0.414		0.856	

Figure 6 shows the overlay of chromatograms obtained from injection of three different brands of syrup samples. No peak of EG and DEG was found in any of these samples. Tables 8 and 9 show the recovery and repeatability (in terms of %RSD) for one of the syrup samples spiked with 200 and 500 µg/mL of EG and DEG, respectively. Excellent recovery and repeatability were achieved for this experiment with syrup sample.

Conclusion

An accurate and rugged method was developed for analysis of ethylene glycol and diethylene glycol in propylene glycol, glycerin, and syrup samples. The LOQ of the method is demonstrated at the 200 µg/mL level in samples. Repeatable results were found for replicates of spiked samples at 200 and 500 µg/mL levels. Good recovery was found at 200 and 500 µg/mL spiked concentration of EG and DEG in different samples. Thus, the method demonstrated in this study proves its usefulness for the routine analysis of excipients and syrup samples for EG and DEG content under the established regulatory limits mentioned in the USP monograph for glycerin³.

References

- 1) Baffi, P.; Elneser, S.; Baffi, M.; De Melin, M. Quantitative Determination of Diethylene Glycol Contamination in Pharmaceutical Products. *J. AOAC Int.* 2000, 83 (4), 793–801.
- 2) U.S. Food & Drug Administration. Lab Method for Gas Chromatography-Mass Spectrometry (GC-MS) Screening Procedure for the Presence of Diethylene Glycol and Ethylene Glycol in Toothpaste. <https://www.fda.gov/food/laboratory-methods-food/lab-method-gas-chromatography-mass-spectrometry-gc-ms-screening-procedure-presence-diethylene-glycol> (accessed 2023-12-01).
- 3) U.S. Department of Health and Human Services, Food and Drug Administration, Center for Drug Evaluation and Research (CDER). Testing of Glycerin, Propylene Glycol, Maltitol Solution, Hydrogenated Starch Hydrolysate, Sorbitol Solution, and other High-Risk Drug Components for Diethylene Glycol and Ethylene Glycol: Guidance for Industry. <https://www.fda.gov/media/167974/download> (accessed 2023-12-01).
- 4) Pharmacopeia. cn USP Method for Glycerol Analysis. http://www.pharmacopeia.cn/v29240/usp29nf24s0_m35420.html#:~:text=Add%2010%20mL%20of%20a,blank%2C%20using%20a%20pH%20meter (accessed 2023-12-01).
- 5) Eccles, R. What is the Role of Over 100 Excipients in Over the Counter (OTC) Cough Medicines? *Lung.* 2020, 198 (5), 727–734. DOI: 10.1007/s00408-020-00390-x.

www.agilent.com/chem

DE87274344

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024
Published in the India, February, 2024
5994-7159EN

