

EVALUATION OF THE D-100 HPLC SYSTEM FOR THE DETERMINATION OF GLYCATED HEMOGLOBIN

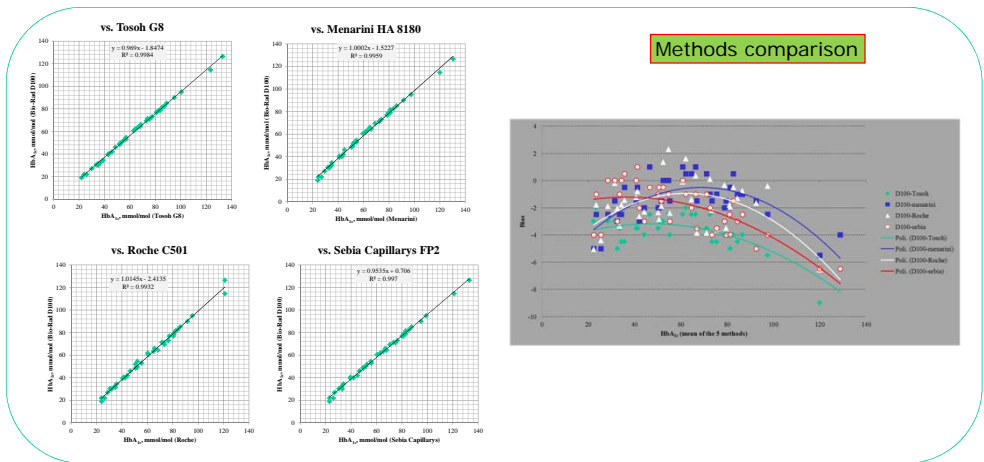
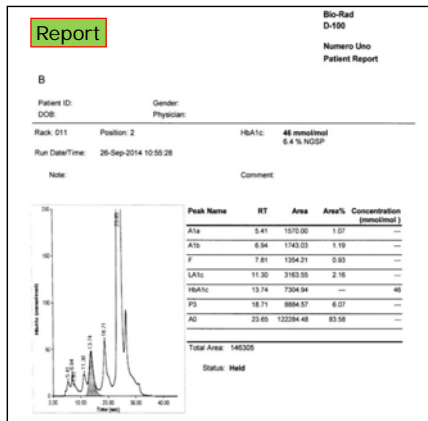
Paleari R¹, Saibene D², Moia R³, Besozzi M³, Varini M⁴, Pezzo M⁴, Ceriotti F², Mosca A¹

¹ Centro per la Riferibilità Metrologica in Medicina di Laboratorio (CIRME), Dipartimento di Fisiopatologia Medico-Chirurgica e dei Trapianti, Università degli Studi, Milano (MI)
² Laboratorio di Standardizzazione, Servizio Medicina di Laboratorio, Ospedale San Raffaele, Milano (MI)
³ Laboratorio di Analisi Cliniche, IRCCS Istituto Auxologico Italiano, Milano (MI)
⁴ Laboratorio Analisi, Ospedale di Suzzara, Suzzara (MN)

Background. The HbA_{1c} testing is set to assume a greater role in the next few years as a consequence of the introduction of HbA_{1c} for the diagnosis in addition to its conventional use for monitoring of diabetic patients, and of the global increasing frequency of diabetes. Accordingly, the methods used for HbA_{1c} determination need to provide excellent performance in terms of analytical quality as well as robustness, usability and throughput. The analytical performance of a new HPLC analyzer, the D-100 System from Bio-Rad Laboratories, has been evaluated

Methods. Precision was tested by using the CLSI-EP5 protocol and measured for 20 working days, 2 runs per day, run in duplicate, on aliquots of frozen blood samples at four different HbA_{1c} levels (30, 47, 62, and 108 mmol/mol). For method comparison, 40 blood samples with HbA_{1c} values well distributed over the measuring interval were analyzed in duplicate according to EP9-A21R. The following comparison methods were used: Menarini HA 8180, Roche Cobas 501, Sebia Capillarys FP2, Tosoh G8. Trueness was evaluated by analyzing two IFCC value-assigned samples. The influence of two common hemoglobin variants (HbS and HbC) was also evaluated.

Results. Total reproducibility was found to be very good at any HbA_{1c} level tested, with CV values always <2 %, well below the recommended goal for imprecision of CV<3%. Method comparison study proved D-100 results were well correlated with those obtained with other methods and provided the following linear regression (least square) equations (D-100 was considered as y, other methods as x), $y=1.000x-1.52$, $R^2=0.995$ (D-100 vs Menarini); $y=1.014x-2.41$, $R^2=0.993$ (D-100 vs Roche); $y=0.953x+0.71$, $R^2=0.997$ (D-100 vs Sebia); $y=0.969x-1.85$, $R^2=0.998$ (D-100 vs Tosoh). With regard to trueness, D-100 presented a bias of -1.2 mmol/mol at HbA_{1c} level of 31.7 mmol/mol and +1.5 mmol/mol at 78.0 mmol/mol respect to the IFCC target values. HbC and HbS were clearly eluted after HbA₀ and not integrated for the calculation of HbA_{1c}.



Precision

Lyophilized samples (Lyphocheck)

HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
35.4	35.4	33.6	31.0	35.6	33.1
Repeatability	0.9	0.7	4.3	2.6	1.9
Between run	0.9	0.7	0.0	0.0	0.6
Between day	0.5	1.1	1.5	1.3	1.2
Total	1.4	1.5	4.5	2.9	2.3

HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
84.1	84.1	79.5	84.7	83.1	81.8
Repeatability	0.4	0.4	1.3	1.4	1.2
Between run	0.0	0.3	0.2	0.8	0.0
Between day	0.2	0.8	1.3	0.8	0.6
Total	0.4	1.0	1.7	1.8	1.3

Fresh human blood

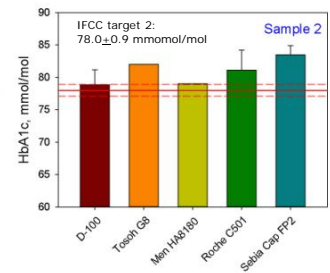
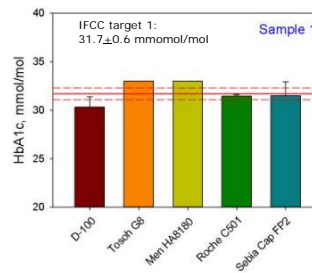
HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
30	30	30	30	30	30
Repeatability	1.0	0.8	2.1	2.1	1.9
Between run	0.5	0.5	0.0	0.0	0.0
Between day	0.4	1.2	0.5	1.7	0.5
Total	1.2	1.5	2.2	2.7	1.9

HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
47	47	47	47	47	47
Repeatability	0.4	0.6	1.9	1.8	1.8
Between run	0.4	0.8	0.0	0.7	0.0
Between day	0.2	0.5	1.2	1.0	0.7
Total	0.6	1.1	2.3	2.2	1.9

HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
62	62	62	62	62	62
Repeatability	0.6	0.5	1.6	1.7	1.2
Between run	0.0	0.7	0.0	0.0	0.0
Between day	0.1	0.7	0.5	0.6	0.7
Total	0.6	1.2	1.6	1.8	1.4

HbA _{1c} mmol/mol	Tosoh G8	Menarini HA 8180	Sebia Cap FP2	Roche C501	Bio-Rad D-100
108	108	108	108	108	108
Repeatability	0.4	0.5	1.1	1.7	1.1
Between run	0.2	0.3	0.0	0.0	0.6
Between day	0.4	0.7	0.6	1.6	0.6
Total	0.5	0.9	1.2	2.3	1.3

Trueness



Analytical interferences

Hb Variant (heterozygosis)	Sample	HbA _{1c} mmol/mol			
		Bio-Rad D 100	Menarini HA8180	Sebia Capillarys	Tosoh G8
Hb C	1	34	34	35	37
	2	39	39	46	41
	3	75	75	89	78
	4	79	79	88	83
Hb D	1	91	65	108	102
Hb S	1	41	38	39	42
	2	42	37	40	43
	3	55	50	55	54
	4	77	76	81	79
(Hb F)	1 (1.0%)	29	30	29	33
	2 (3.5%)	35	36	32	39
	3 (1.0%)	36	34	nd	36
	4 (2.2%)	41	36	nd	39
	5 (2.6%)	40	35	nd	39

Method	HbA _{1c} mmol/mol			
	Influence of triglycerides		Influence of bilirubin	
	Basal	Spiked	Basal	Spiked
Bio-Rad D100	40	40	42	43
Menarini HA8180	40	40	41	41
Sebia Capillarys 2FP	39	38	39	38
Tosoh G8	43	43	45	45

Conclusions. The Bio-Rad D-100 system is a fully automated, user-friendly, high throughput HPLC system giving accurate and reproducible results. The system showed a reduced hands on time, a simple calibration process, a very good workflow efficiency thanks to the intuitive interface, and an high speed HbA_{1c} assay in 45 sec. Bio-Rad D-100 therefore displays the appropriate characteristics to be used as a routine method in clinical laboratories.