

Diabetes Correlation Card

NGSP A1C% ¹	eAG (mg/dL) ²
4.0	68
4.5	82
5.0	97
5.1	100
5.2	103
5.3	105
5.4	108
5.5	111
5.6	114
5.7	117
5.8	120
5.9	123
6.0	126
6.1	128
6.2	131
6.3	134
6.4	137
6.5	140
6.6	143
6.7	146
6.8	148
6.9	151
7.0	154
7.1	157
7.2	160
7.3	163
7.4	166
7.5	169
7.6	171
7.7	174
7.8	177
7.9	180

(See back for references)

diabetes.bio-rad.com



Diabetes Correlation Card



NGSP A1C% ¹	eAG (mg/dL) ²
8.0	183
8.1	186
8.2	189
8.3	192
8.4	194
8.5	197
8.6	200
8.7	203
8.8	206
8.9	209
9.0	212
9.1	214
9.2	217
9.3	220
9.4	223
9.5	226
9.6	229
9.7	232
9.8	235
9.9	237
10.0	240
10.1	243
10.2	246
10.3	249
10.4	252
10.5	255
10.6	258
10.7	260
10.8	263
10.9	266
11.0	269
11.5	283
12.0	298
12.5	312

^{1,2} American Diabetes Association
 <<http://professional.diabetes.org/glucosecalculator.aspx>>

Bio-Rad A1c • Be the difference