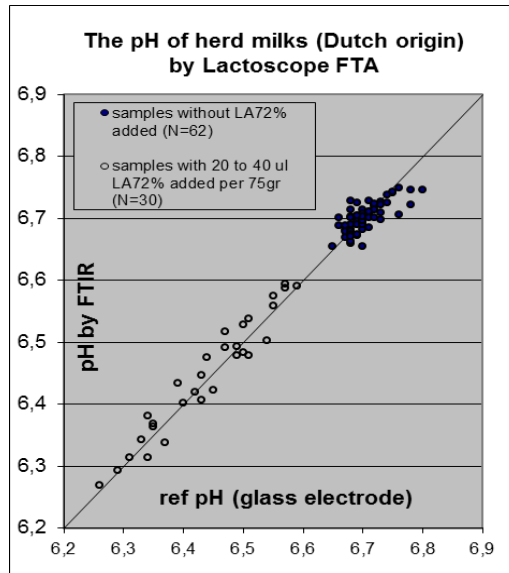
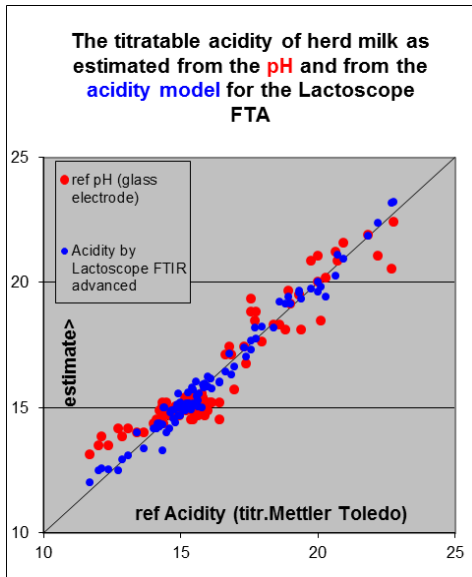
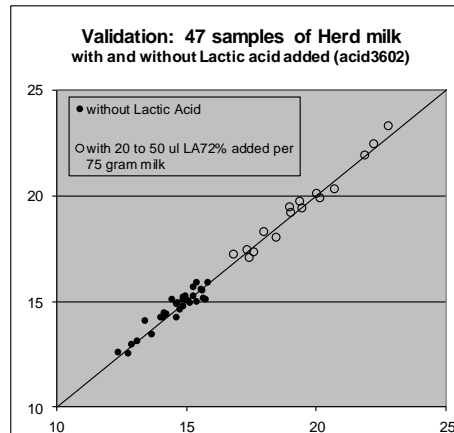




Author: WB		<b>Type: application sheet</b>				
Date	27-05-2008	class	VL	Mod.code	Units	instrument
Rev.	acidity1.0	type	PLS	Parm 1501	Titrateable Acidity of milk in mmol lactic acid/Liter	FTA
Doc. contr	ensur	Comm.	public	Parm 151		Combi FTIR 400
		Validation notes	0			



ref Acidity from:	N=92		
Prediction function	<b>SEC</b>	<b>CV</b>	<b>R<sup>2</sup></b>
<b>Y= a ref pH+b</b>	0,78	4,8%	0,90
<b>Y= a * Acid IR</b>	0,35	2,1%	0,99



N	Mean	CV	SEP
47	16,2	2,0%	0,33

