

---

# **IgG core a-fucosylation and its impact on Fc $\gamma$ RIIIa binding**

**MipTec 21.09.2011**

***Sandra Grau***

***Roche Glycart AG***

---

# Introduction

---

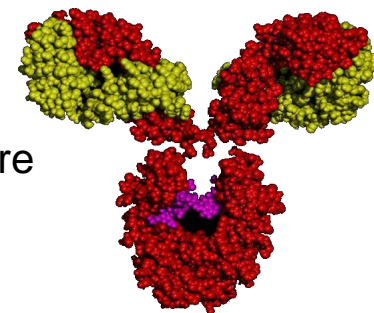
## TagLite Results

## Outlook

# Introduction

## *Antibody therapies*

Roche



- Monoclonal antibodies represent a growing class of therapeutics with more than 20 molecules licenced for the treatment of cancer and chronic diseases.
- Major indications: oncology, infectious diseases and autoimmunity.
- Efficacy results from their specificity to the antigen target as well as the activation of effector functions.
- In oncology one relevant mechanism of action is antibody-mediated cellular cytotoxicity (ADCC):
  - This was shown for various antibody therapeutics such as rituximab (anti-CD20) and trastuzumab (anti-Her2).
  - The mechanism underlying ADCC is the binding of FcγRIIIa on natural killer cells (NK cells) to the Fc portion of an IgG and killing of the target cell

# Introduction

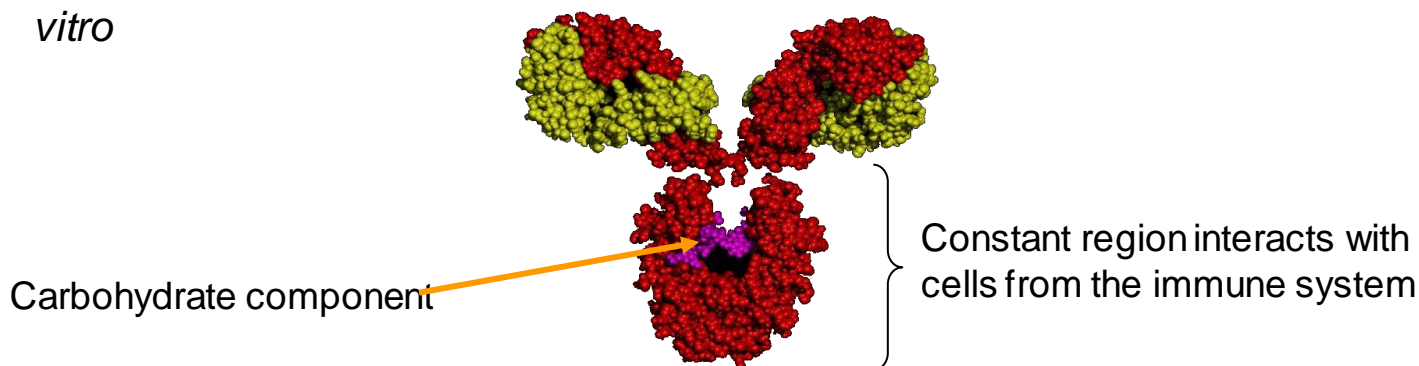
## Glycoengineering

**Main goal for next generation therapeutic antibodies was:  
Increase the binding of the antibody to activating  $\text{Fc}\gamma\text{Rs}$  ( $\text{Fc}\gamma\text{RIIIa}$ ).**

➤ *Two different strategies:*

- Amino acid mutations in the Fc part of the antibody
- Changing the carbohydrate moieties in the Fc portion of the antibody

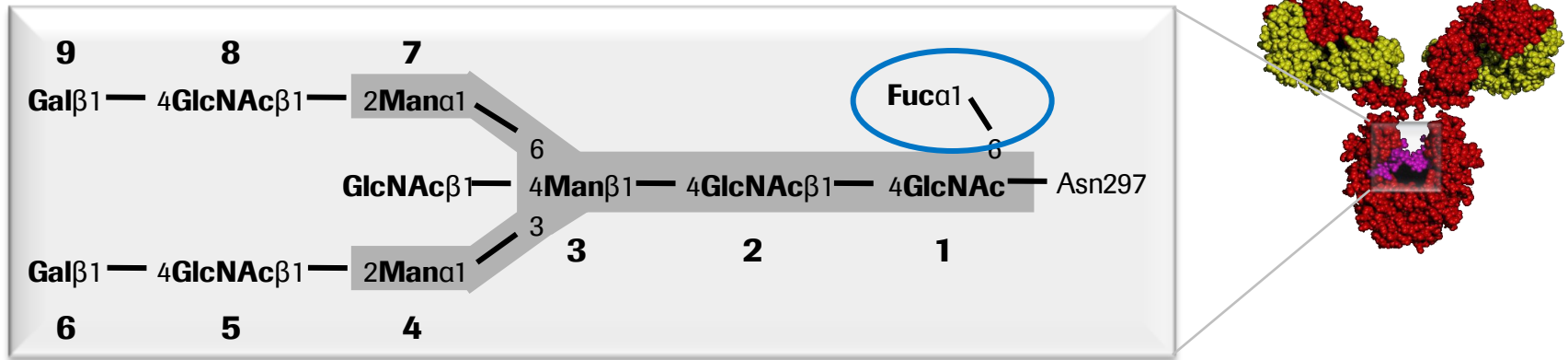
→ a-fucosylated antibodies show increased binding to  $\text{Fc}\gamma\text{RIIIa}$  and enhanced ADCC *in vitro*



# Introduction

## Glycoengineering

### ➤ Glycosylation of IgGs



**Removal of fucose → increased affinity for FcγRIIIa**

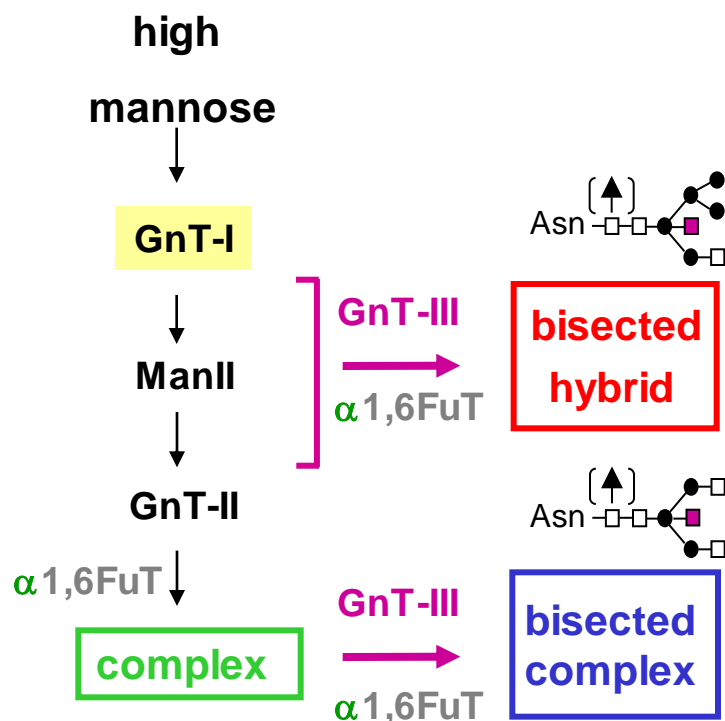
**How can one generate antibodies  
lacking the fucose?**

# Introduction

## Glycoengineering

### → Engineering cell lines overexpressing GnTIII

#### N-linked glycosylation



<b>GnT-III</b>	→	<b>bisected</b>
competes against		
$\alpha 1,6$ -FuT	→	non-fucosylated
<b>ManII</b> and <b>GnT-II</b>	→	<b>hybrid</b>

**Why has the presence of the fucose such a great impact on binding?**

□ N-acetylglucosamine (GlcNAc)

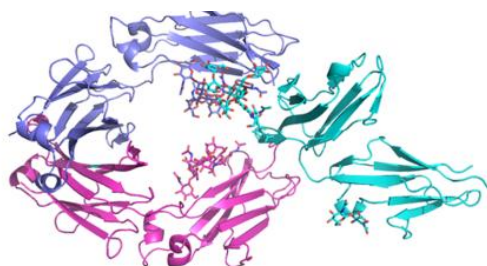
▲ Fucose

● Mannose

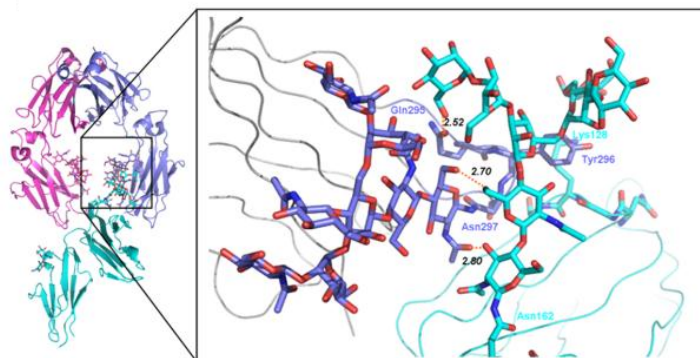
# Introduction

## *Crystal structure of Fc $\gamma$ RIIIa complexed with either $\alpha$ -fucosylated or fucosylated Fc*

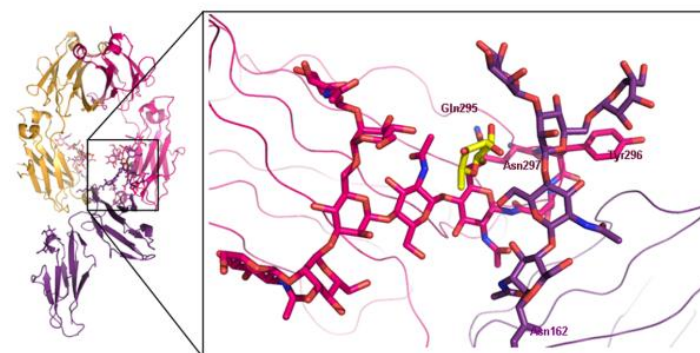
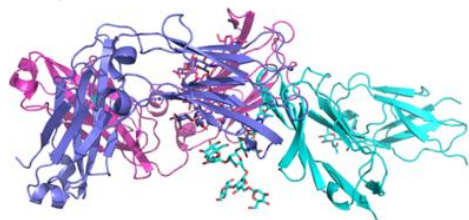
***Fc $\gamma$ RIIIa /  $\alpha$ -fuc Fc***



CH3 CH2 hFc $\gamma$ RIIIa



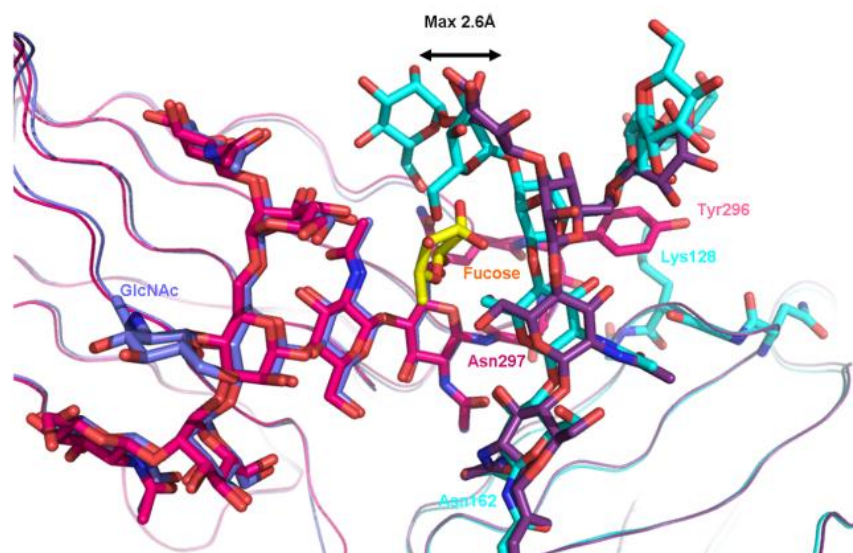
***Fc $\gamma$ RIIIa / fuc Fc***



- Carbohydrate-carbohydrate mediated interactions are responsible for an up to 100-fold gain in binding affinity for  $\alpha$ -fucosylated vs. fucosylated IgGs.

# Introduction

## ***Crystal structure of Fc $\gamma$ RIIIa complexed with either $\alpha$ -fucosylated or fucosylated Fc***



- The Fc-core fucose (highlighted in yellow) has to accommodate in the interface and the Asn162-receptor glycan has to move ( $\leftrightarrow$ ).
- The result is a direct, steric inhibition caused by core fucose for the carbohydrate-mediated interaction with Fc $\gamma$ RIIIa.
- The structures provide a molecular mechanism explaining the increased affinity for the receptor of  $\alpha$ -fucosylated antibodies.



# Introduction

## *Characterization of monoclonal IgGs*

### Carbohydrate analysis

High throughput  
ProtA purification



Endoglycosidase  
digest



MALDI TOF MS  
analysis

### Biological activity

ADCC

### Affinities

Surface Plasmon  
Resonance

**Could the TagLite technology combine  
some of these analysis?**

# Introduction

---

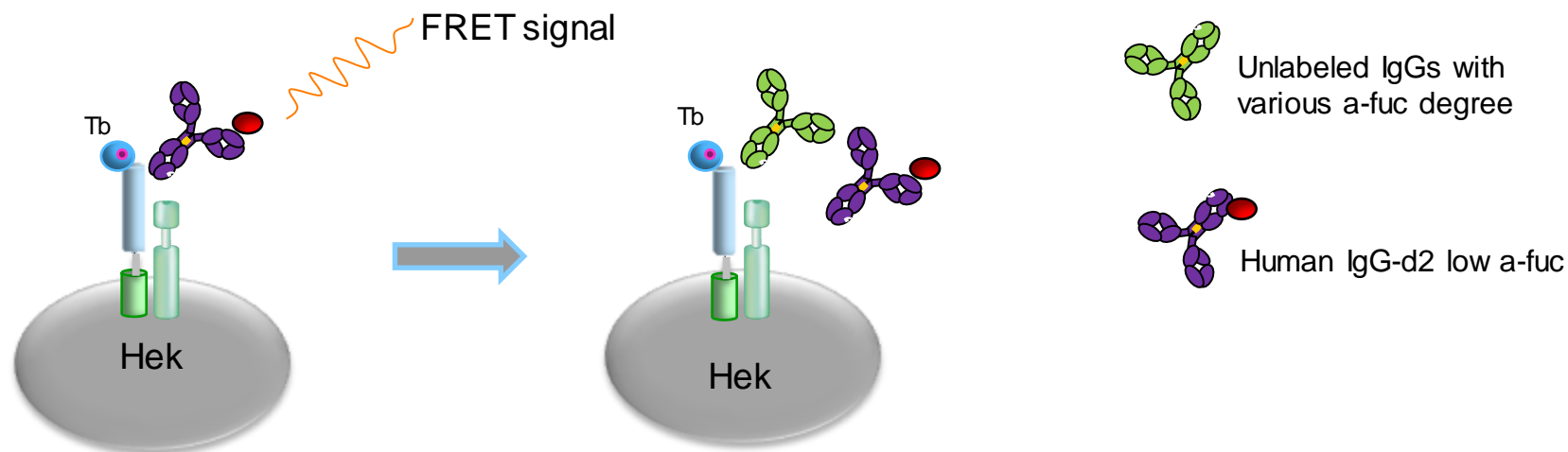
## TagLite Results

---

## Outlook

# TagLite – results

## *Fc $\gamma$ RIIIaV158 and F158 competition assay*

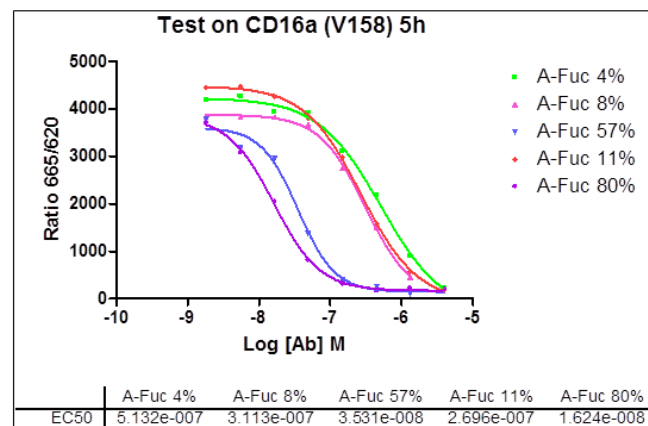
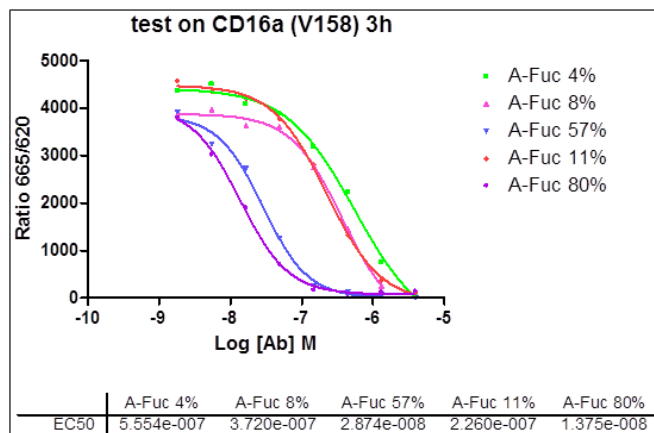
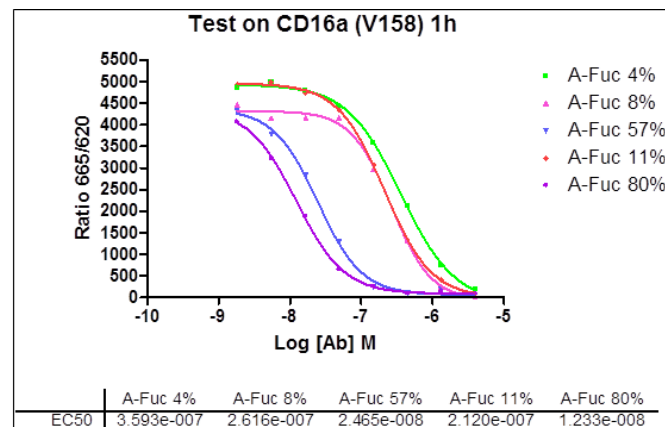
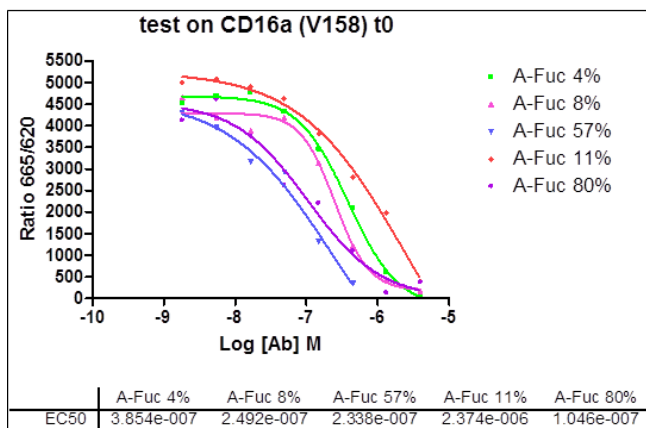


### **Assay setup:**

- 10.000 cells per well (expressing huFc $\gamma$ RIIIaV158 or F158 labeled with Tb)
- IgGs with various a-fucosylation degrees (4, 8, 11, 57 and 80 %) conc. 4000 to 1.8 nM final in well (1:3 dilutions).
- Labeled IgGs (human IgG-d2); conc. 50 nM final in well
- Incubation time: 0h, 1h, 3h, 5h @ RT

# TagLite – results

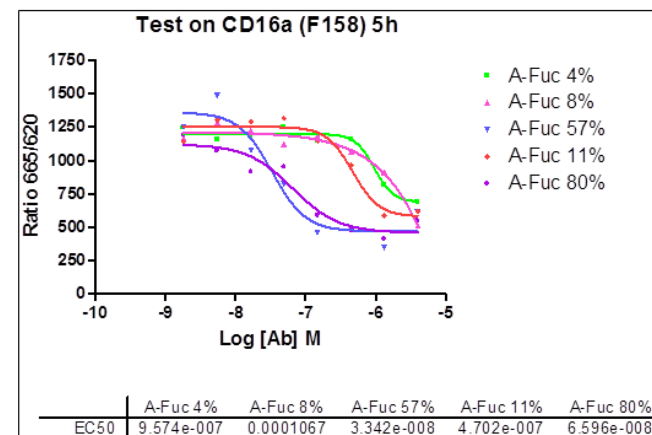
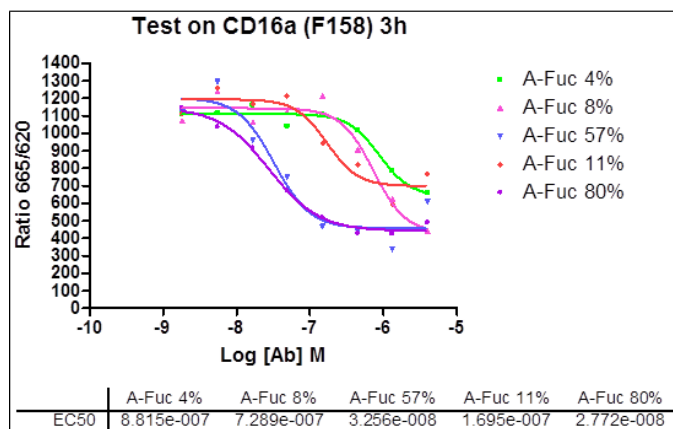
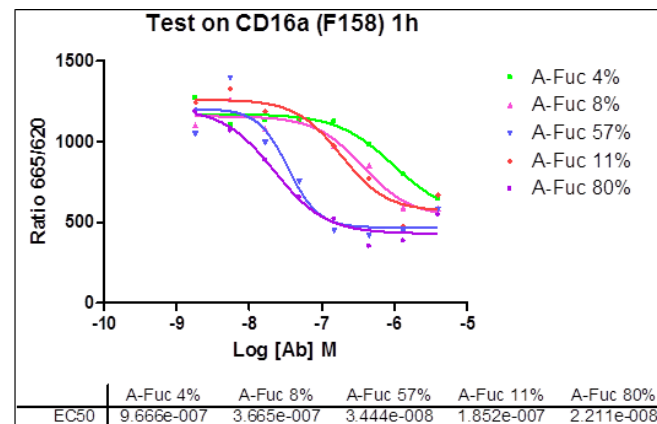
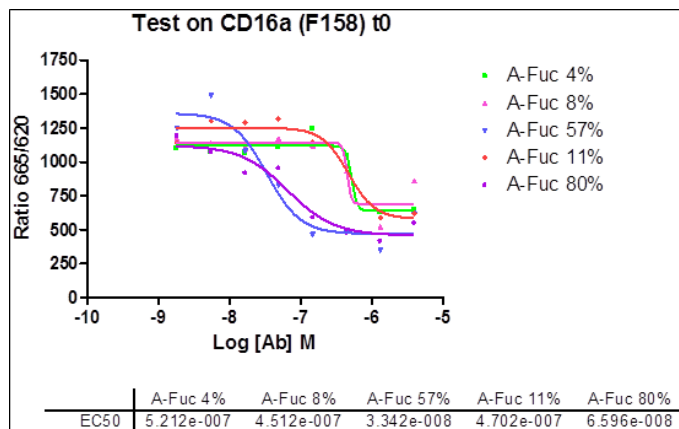
## *Fc $\gamma$ R11a(V158)*



- The assay works nicely with the high affinity receptor.
- Already after one hour incubation ranking of the abs can be detected.
- The 2 IgGs with high a-fuc degree compete much better than the 3 IgGs with lower a-fuc degree.

# TagLite – results

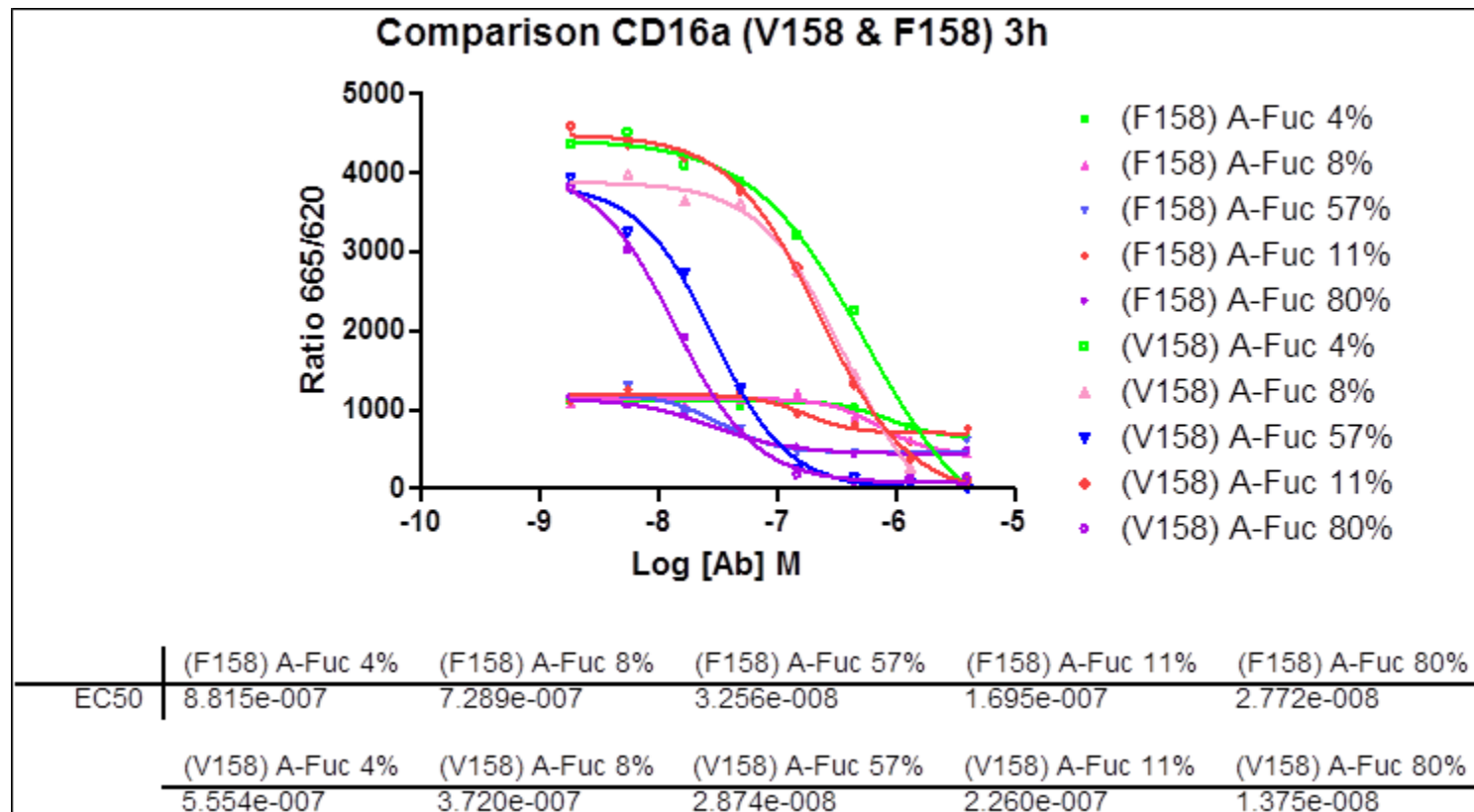
## *Fc $\gamma$ R11a(F158)*



- The assay also works with the low affinity receptor but the assay window is much smaller.
- Still also here the higher a-fucosylated IgGs compete better than the ones with lower a-fuc content.

# TagLite – results

## Comparison $Fc\gamma RIIIa(V158)$ and $Fc\gamma RIIIa(F158)$



# TagLite – results

## Correlation with other methods



**huFcγRIIIaF158**

	MALDI	Biacore	TagLite
	a-fuc (%)	KD (nM)	EC50 (nM)
832	4	1300	881
834	8	940	728
835	11	845	170
836	57	68	32
842	80	39	27

a-fuc increase ↓ improve in KD/EC50 ↓

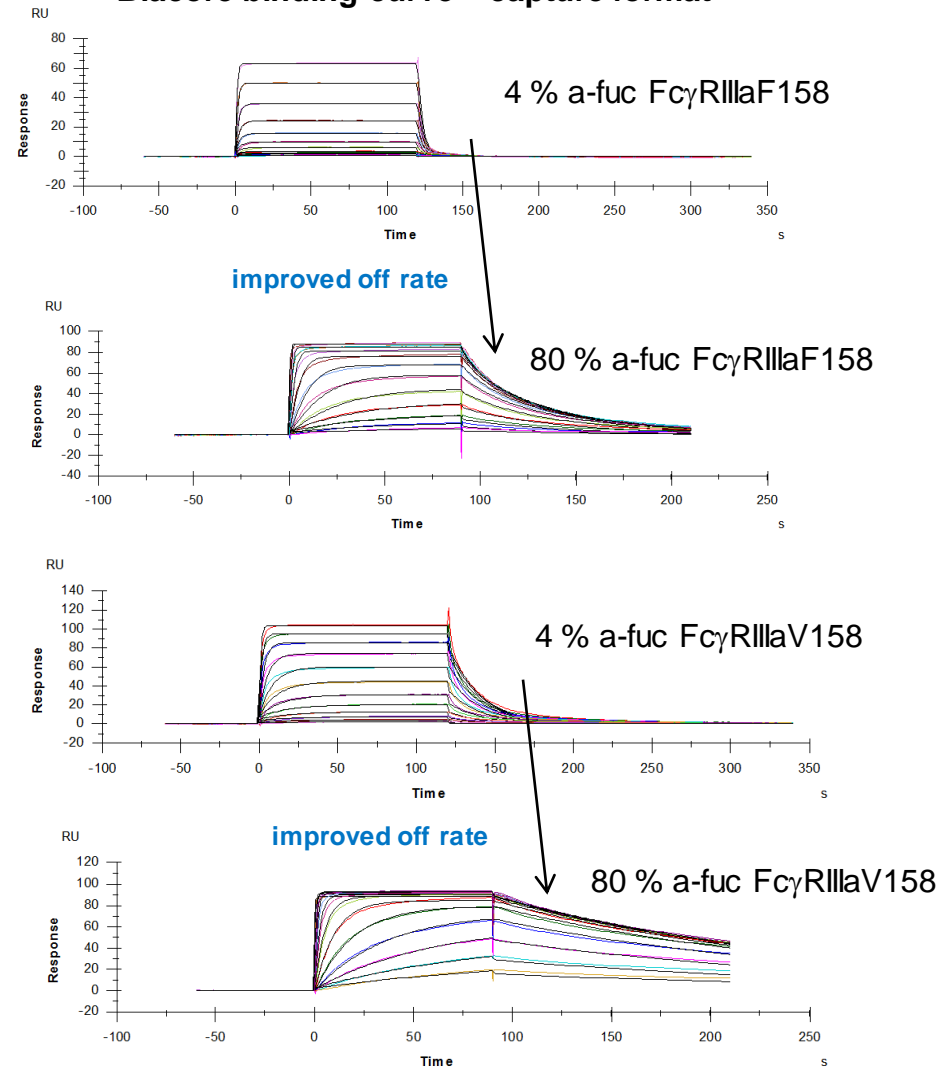
**huFcγRIIIaV158**

	MALDI	Biacore	TagLite
	a-fuc (%)	KD (nM)	EC50 (nM)
832	4	200	555
834	8	201	370
835	11	111	226
836	57	17	28
842	80	7	13

a-fuc increase ↓ improve in KD/EC50 ↓

Note: KD values (Biacore) depend a lot on the format used.  
In the present format KD values for a-fuc abs tend to be lower than with other formats. You can expect values ranging from 500-700 nM for FcγRIIIaV158 and low a-fuc abs.

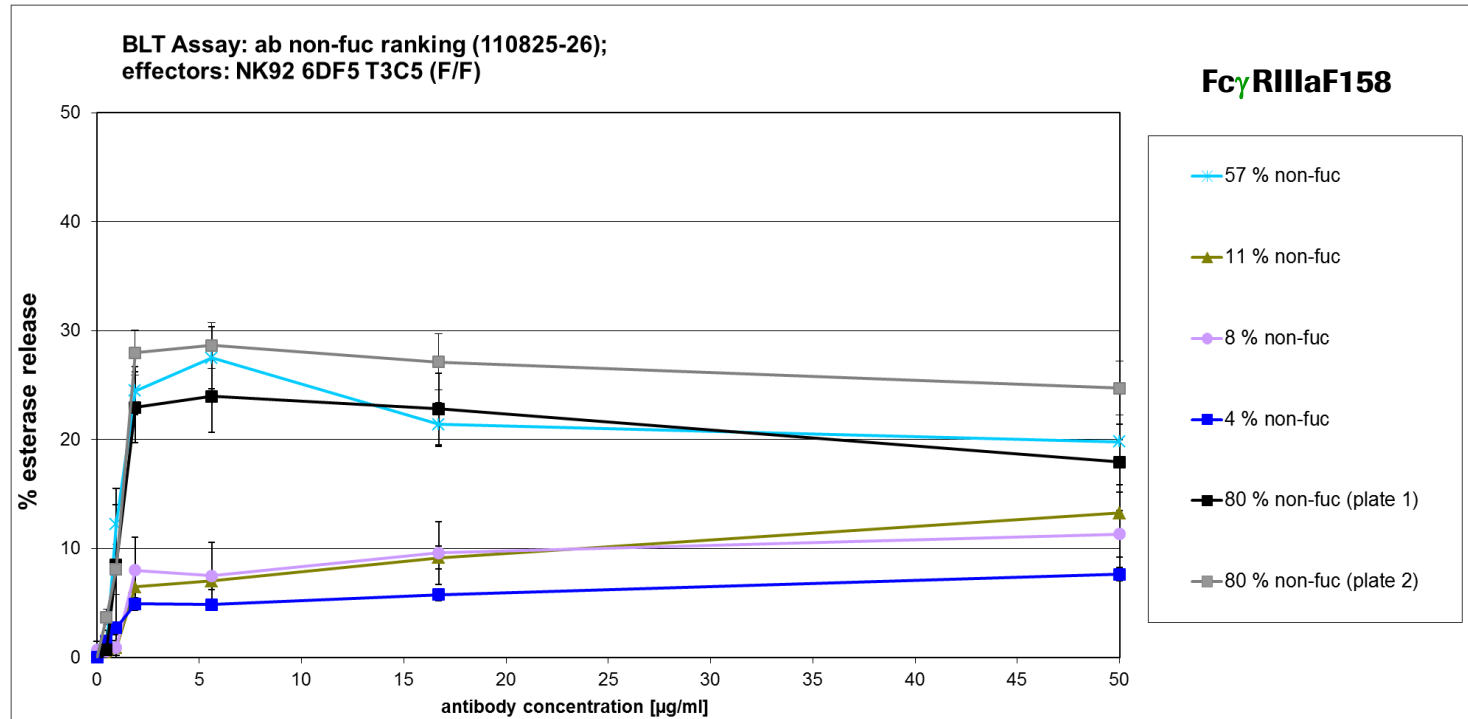
Biacore binding curve – capture format



# TagLite – results

## *Correlation with other methods*

### Biological activity - BLT assay



- The BLT assay measures the release of esterase upon activation of NK cells
- The graph shows that the 2 IgGs containing higher a-fuc degree lead to a 3 fold higher release of esterase than the ones with lower a-fuc degree.
- The release correlates well with killing of the target cell.



# Introduction

## TagLite Results

---

## Outlook

---

- The TagLite technology provides a nice, robust and easy to use tool to study binding of antibody Fc portions to Fc $\gamma$ RIIIa on cells.
- The data correlate well with data from Biacore and cell based assays like ADCC
- An advantage compared to Biacore analysis is that binding occurs with native receptor embedded in the cell membrane rather than purified soluble forms which translates better into cell based assays.
- Interesting would be to test also other hu Fc $\gamma$  receptors as well as muFc $\gamma$ R which might be of importance for mouse models.
- Also to be checked: Could you determine a-fucosylation degree simply by using a standard curve of IgGs with various a-fucosylation degree and therefore use the TagLite assay for screening purposes?

# Acknowledgements

## **Roche Glycart**

Pablo Umana

Christiane Jaeger

Claudia Ferrara

Sylvia Herter

## **Cisbio Bioassay**

Delphine Jaga

Hamed Mokrane



*We Innovate Healthcare*