



## Targeting gene dependencies in MYC overexpressing Multiple Myeloma

Lama Hasan Bou Issa<sup>\* 1</sup>, Léa Fléchon<sup>1</sup>, William Laine<sup>1</sup>, Aicha Ouelkdite<sup>1</sup>, Guillaume Escure<sup>2</sup>, Silvia Gaggero<sup>1</sup>, Tiziano Ingegnere<sup>1</sup>, Romanos Sklavenitis Pistofidis<sup>3</sup>, Irene M. Ghobrial<sup>3</sup>, Thierry Facon<sup>2</sup>, Suman Mitra<sup>1</sup>, Jérôme Kluza<sup>1</sup>, Bruno Quesnel<sup>1, 2</sup>, Salomon Manier<sup>1, 2</sup>

<sup>1</sup>Univ Lille, Canther, INSERM UMR-S1277 - CNRS UMR9020, <sup>2</sup>CHU Lille, Department of Hematology, Lille, France,

<sup>3</sup>Dana-Farber Cancer Institute, Harvard Medical School, Boston, MA, United States of America

**Background.** Multiple myeloma (MM) is an incurable plasma cell malignancy in which *MYC* alterations play an essential role in the progression and correlate with adverse outcome. However, *MYC* is not therapeutically targetable.

**Hypothesis.** To overcome this, we hypothesized that the proliferative advantage promoted by *MYC* overexpression induces differential genomic dependencies that can be exploited therapeutically.

**Methods.** We searched for genomic vulnerabilities associated with *MYC* overexpression using genome-scale pooled short hairpin RNA screening data (Project Achilles). We generated an isogenic model of *MYC* overexpression (OE) in U266 cell line. We performed RNA-seq, quantitative proteomics and a drug screening of ~2000 small molecules. For validation, we performed pharmacological inhibition of glutaminolysis and shRNA-mediated *GLS1* knockdown. CellTiter-Glo® was used to assess viability in both drug screen and validation tests. To understand the differential metabolic rewiring, we performed metabolomic and Seahorse XF analysis.

**Results.** Achilles analysis revealed main dependencies associated with *MYC* overexpression specifically *GLS1* (glutaminase) and *SLC1A1* (glutamine transporter). Our drug screen spotlighted NAD synthesis inhibitors and mTORC1, which rely on intracellular glutamine pool, had preferential effect on U266/*MYC*. Both RNA-seq and quantitative proteomics showed no significant upregulation of glutaminolysis-related genes, suggesting a non-oncogenic dependency. Our data showed that *MYC* OE cells failed to proliferate under glutamine deprivation and had higher sensitivity to CB-839 and V-9302 inhibiting *GLS1* and *SLC1A5*, respectively. Using the Seahorse Bioanalyzer, we observed higher oxygen consumption rate (OCR) in U266/*MYC* which was induced by glutamine. Furthermore, CB-839 had a potent effect on U266/*MYC* observed at basal and maximal OCR highlighting the glutamine role controlling mitochondrial OXPHOS in *MYC* OE cells. The metabolomic analysis showed higher *GLS1* activity in U266/*MYC* with elevated glutamine to glutamate flux. Notably, CB-839 in U266/*MYC* results in more pronounced changes in TCA cycle and energy debt. This effect was blunted by co-incubation with  $\alpha$ KG. Interestingly *GLS1* inhibition was not limited to this, but extended to redox balance and amino acid biosynthesis.

**Conclusion.** A comprehensive understanding of these vulnerabilities can provide a powerful roadmap to guide innovative therapeutic approaches. In this prospective, our results are important in finding new strategies to target *MYC* in clinic.